



Directions for Use

Standard Filter Cartridges and Mini Cartridges

1. Labeling

Product name, pore size or retention rate, lot number and order code are indicated on the top adapter of the standard filter cartridges.

The label on the plastic bag and the outer packaging includes the product name, pore size or retention rate, expiry date and lot number along with the order code.

2. Installing the Filter Cartridges

- Cut open the plastic bag at the bottom end of the filter cartridge (on the O-ring side).

⚠ Important

Wet the gaskets with water or the medium to be filtered!

- Leave the cut-open filter bag as a protection around the filter cartridge while installing it.

⚠ Important

During installation, avoid any damage to the adapter or O-ring! For this purpose, only touch the filter cartridge in the bottom area.

- Insert the filter cartridge into the bottom of the housing using a slight rotating movement.
- Remove the protective bag.
- Close the housing by installing the bell on the top of the base.

⚠ Important

If you cannot use standard silicon O-rings, replace them with equivalent O-rings before installing the filter cartridge.

3. Wetting and Rinsing (in the Direction of Filtration)

⚠ Important

Perform a rinsing cycle before every sterilization routine and integrity testing (membrane filters only). This will help to ensure a filtration system that is as particulate-free as possible after installing the cartridges.

This step is not necessary for hydrophobic filter elements (e.g. Sartofluor®).

- For this purpose, open the vent valve during filling.
- Fill the cartridge housing with de-ionized water, filter medium or, for critical applications, with pyrogen-free distilled water, until liquid flows out of the vent valve.

⚠ Notice

If the filter cartridge is a product with a PTFE membrane (e.g. Sartofluor® GA, Sartofluor® LG or Aerosart), it has to be wetted using a liquid that has surface tension < 28 dyne/cm.

⚠ Important

By keeping the vent valve slightly open during the first minute of the rinse cycle, you ensure that the air is removed and the filter is properly wetted. No air remains in the housing.

- Close the vent valve.
- Rinse at a differential pressure of 0.3 bar for approx. 3–5 minutes.
- After rinsing, open the vent and the run-off drain valve.
- Drain the cartridge housing. This can also be done with compressed air or nitrogen at a maximum inlet pressure of 0.5 bar with the vent valve closed.
- The cartridge is properly wetted once you have carried out the rinsing procedure as described above.

4. Sterilization

Autoclaving

Sterilization temperature: Up to 134°C, 2 bar

- Autoclave the filter cartridges for a minimum of 30 minutes after the sterilization temperature has been reached. Cover inlet and outlet and any opened valves on the housing with autoclave paper to form a vapor permeable seal or aluminum foil.

In-Line Steam Sterilization

Saturated Steam: Up to 134°C, 2 bar

- Open the vent valve, run-off valve and, if appropriate, the check valve.
- Let the steam flow in slowly and monitor that the differential pressure of 0.5 bar is not exceeded during the pressurization phase.
- Close almost all valves as soon as steam escapes from them.
- Set the desired sterilization steam pressure at the filter outlet through the different valves (steam, vent, check and outlet valves).

⚠ Important

The differential pressure between the inlet and outlet sides should not exceed 0.5 bar; this will ensure the longest service life possible!

- Steam sterilize the filter cartridges (once the sterilization temperature has been reached at the outlet) in or against the filtration direction for 30 minutes under at the most constant pressure & temperature conditions possible.
- After completing the sterilization cycle, turn off the steam supply.
- Wait until no more steam escapes from the outlet and the inlet and outlet pressures have dropped to atmospheric.
- Close the valves on the sterile side.

⚠ Notice

During the cooling phase, avoid a vacuum condensation as follows: After steam sterilization, cold-blow the filter system. For this purpose, use a supporting gas pressure (compressed air, carbon dioxide or nitrogen). The pressure must be 200 mbar higher than the steam pressure applied during sterilization.

- With closed upstream housing valves, you can now carry out an integrity test.

Sanitization with Hot Water (Food & Beverage Application)

In selected applications, sanitization can be carried out by circulating hot water through the system. Make sure that the housings are filled completely.

- During filling, leave the vent valve open.
- The sanitization time takes 30 minutes after the water at the outlet has reached a minimum of 85°C.
- After completing sterilization, turn off the hot water supply.

⚠ Notice

During the cooling phase, avoid a vacuum condensation as follows: After sanitization, cold-blow the filter system. For this purpose, use a supporting pressure (compressed air, carbon dioxide or nitrogen) that is 200 mbar higher than the atmosphere pressure applied during sterilization.

- With the upstream housing valves closed, you can now carry out an integrity test.

5. Integrity Tests and Integrity Test Data (only membrane filter cartridges)

After wetting and cooling the membrane filter cartridges, the integrity can be tested.

- Carry out this test with an automatic integrity testing device such as Sartocheck® or using a manual integrity.

For more information about integrity testing and product-specific integrity test data, please refer to the “Product Data” file on the enclosed CD.

6. Maximum Allowable Differential Pressures

In the filtration direction:

- At 20°C max. 5 bar
- At 80°C max. 2 bar

Against the filtration direction:

- At 20°C max. 2 bar

7. Filtration

To make use of the entire filter surface, vent the housing completely at the start of filtration. For this purpose, briefly open the vent valve until the liquid to be filtered escapes and then set the desired filtration pressure.

8. Changing the Filter Cartridges

At the latest, replace the filter cartridges as part of the liquid filtration procedures whenever

- the maximum allowable differential pressure has been reached
- or
- the flow rate drops significantly.

9. Applications Support

If needed, please request the corresponding documents with further technical data, on how to carry out product integrity tests or special information about additional applications. Moreover, if you have specific questions about product validation, you can also contact our Validation Service or our Technical Support departments; they will be happy to assist you with all technical questions and optimization procedures.

10. Returning Used Filter Elements

⚠ Important

Before returning used filter elements to Sartorius Stedim Biotech for an additional testing, make sure that they have been sterilized and decontaminated properly. This has to be certified in the required Return Shipment Form that you can request from your local Sartorius Stedim Biotech representative. Otherwise, the provisions of the Employment Protection Act prohibit us from processing the returned products.

11. Liability

Sartorius Stedim Biotech does not assume responsibility in the event of improper use of the filter cartridges. In the interest of further developing Sartorius Stedim Biotech products, we reserve the right to make changes to the product.

⚠ Notice

The enclosed CD contains PDF files in different languages with further product-specific parameters for the filtration products (refer to “CD Information” and “CD Installation” on the next page). Please read this information carefully before using the product.

Directions for Use

Maxicaps[®], Midicaps[®] and Capsule

1. Labeling

Product name, pore size or retention rate, order code, lot number, operating parameters, operating direction and expiry date are indicated on the housing of the Maxicaps[®], Midicaps[®] and the capsules. Product name, pore size or retention rate, order code, lot number and the consecutive individual item number are indicated on an outer sleeve of the Maxicaps[®] banderole.

The operating parameters (max. operating temperature) and the operating direction are also indicated. The label on the plastic bag and the outer packaging includes the product name, pore size, expiry date and batch number along with the order number.

⚠ Notice

Selected products (with "Sterile product" on the label) are heat-sterilized in the primary packaging using a process validated in accordance with DIN/EN 554. Sterility is only ensured if the primary packaging is intact. For that reason, always check to make sure that the packaging is intact before using a sterile product. If the packaging is damaged, the user must sterilize the product as described in Section Autoclaving before using the product as a sterile filter.

2. Wetting and Rinsing (in the Filtration Direction)

⚠ Important

Do not operate venting valves when the differential pressure is higher than 500 mbar.

⚠ Important

Make sure the valves are closed before wetting and rinsing the Maxicaps[®], Midicaps[®], and capsules; this will prevent any liquid from escaping from the valves unintentionally.

⚠ Important

Carry out a rinsing cycle before every sterilization routine and integrity test. This will help to ensure a filtration system, that is as particulate free as possible after the installation.

- For this purpose, open the vent valve during filling.
- Fill the housing with de-ionized water, filter medium or, for critical applications, with pyrogen-free distilled water, until liquid escapes from the vent valve.

⚠ Notice

If the filter device cartridge is a product with a hydrophobic membrane (e.g. Sartofluor[®] or Sartopore[®] Air), it has to be wetted using a liquid with a surface tension < 28 dyne/cm.

⚠ Important

By keeping the vent valve(s) slightly open during the first minute of the rinse cycle, you ensure that no air remains in the housing.

- Close the vent valve.
- Rinse at a differential pressure of 0.3 bar for approx. 3–5 minutes.
- After rinsing, open the vent and the run-off valve.
- Drain the housing. This can also be done with compressed air or nitrogen at a maximum inlet pressure of 0.5 bar with the vent valve closed.
- The filter device cartridge is properly wetted once you have carried out the rinsing procedure as described above.

3. Sterilization

Autoclaving

⚠ Important

Gamma irradiatable filter devices only may be sterilized by autoclaving three times at the most!
For other products refer to product datasheet.

Sterilization temperature: Up to 134°C, 2 bar

- Autoclave the filter device cartridges for a minimum of 30 minutes after the sterilization temperature has been reached. Cover inlet and outlet and any opened valves on the housing with autoclave paper or aluminum foil to form a vapor-permeable seal.

⚠ Notice

Pay attention that, during autoclaving, the adapter parts (e.g. stainless steel parts) connected to the Maxicaps[®], Midicaps[®] or shroud are not overtightened. Also, prevent heavy connecting parts from exerting traction on the device's product housing as this may otherwise cause deformation.

In-Line Steam Sterilization

⚠ Important

Maxicaps[®], Midicaps[®] and Capsules may not be in-line steam sterilized!

CD Information

The enclosed CD contains the operating instructions as a PDF file. If the CD is missing, you can obtain a copy from us by specifying the order number:

Material No. 85032-539-79
Publication No. SPK6168cc

Sartorius Stedim Biotech GmbH
Technical Editorial Department
August-Spindler-Strasse 11
37079 Goettingen, Germany
tech.pubs@sartorius.com
www.sartorius.com

Gamma Sterilization

⚠ Important

Gamma irradiatable filter devices can be sterilized one time by gamma irradiation with max. 50 K Gy. Only one sterilization method (autoclaving or gamma sterilization) has to be performed.

4. Integrity Tests and Integrity Test Data

After wetting and sterilizing, Maxicaps[®], Midicaps[®] and Capsules with a membrane can be integrity tested after cooling down to room temperature.

- It is highly recommended that testing is performed by an automatic integrity testing device like Sartocheck[®].

For more information about integrity testing and product-specific integrity test data, please refer to the "Product data" file on the enclosed CD.

5. General Safety Instructions

When installing Maxicaps[®], Midicaps[®] or Capsules, ensure that the capsule is suitable for its intended purpose and that it is installed correctly. External mechanical impacts caused by improper handling of Maxicaps[®], Midicaps[®] and Capsules on site during or after unpacking, e.g. capsule fallen down or obvious mechanical shock and stress to the capsule housing, must be avoided. Such external mechanical impact can lead to cracked and damaged Maxicaps[®], Midicaps[®] and Capsules. When pressurized during filtration or integrity testing, this can result in leakage or rupture failures. For this reason, Maxicaps[®], Midicaps[®] and Capsules must no longer be used, under any circumstances, once they have been subjected to such external mechanical impacts.

General rules and safety instructions when working with pressurized systems must be followed.

6. Maximum Allowable Differential Pressures

In the filtration direction:

- At 20°C max. 4 bar (Capsules, Maxicaps[®])
- At 20°C max. 5 bar (Midicaps[®], T-Style Maxicaps[®])

Against the filtration direction:

- At 20°C max. 2 bar

7. Filtration

To make use of the entire filter surface, vent the housing completely at the start of filtration. For this purpose, briefly open the vent valve until the liquid to be filtered escapes. Close the vent valves and increase pressure to the desired filtration pressure.

8. Product Change

At the latest, replace the filter cartridges as part of the liquid filtration procedures whenever

- the maximum allowable differential pressure has been reached
- or
- the flow rate drops significantly.

9. Applications Support

If needed, please request the corresponding documents with further technical data, on how to carry out product integrity tests or special information about additional applications. Moreover, if you have specific questions about product validation, you can also contact our Validation Service or our Technical Support departments; they will be happy to assist you with all technical questions and optimization procedures.

10. Returning Used Filter Elements

⚠ Important

Before returning used filter elements to Sartorius Stedim Biotech for additional testing, make sure that they have been sterilized and decontaminated properly. This has to be certified in the required Return Shipment Form that you can order from your local Sartorius Stedim Biotech representative. Otherwise, the provisions of the Employment Protection Act prohibit us from processing the returned products.

11. Liability

Sartorius Stedim Biotech does not assume liability if the filter cartridges are handled improperly. In the interest of further developing Sartorius Stedim Biotech products, we reserve the right to make changes to the product.

⚠ Notice

The enclosed CD contains PDF files in different languages with further product-specific parameters for the filtration products (refer to "CD Information" and "CD Installation" on this page). Please read this information carefully before using the product.

System requirements:

- Windows[®] XP or higher
- Adobe[®] Reader[®] 5.0 or higher

CD Installation

Proceed as follows to use the PDF files:

- Start Windows.
- Insert the CD in the CD drive.
- The software should start automatically.
- If not, run Start.exe on the CD to start the application.

On other operating systems please use start.html. To eject the drive you need to close the browser window (Internet Explorer).

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0
www.sartorius.com

Specifications subject to change without notice.
Copyright Sartorius Stedim Biotech GmbH.
DIR: 2622109-000-00
Status: 08 | 2020

List of Sartorius material numbers applying to EPA-FIFRA

16702-27	5102507H4	5107307H1-SS
16702-28	5102507H4--EP	5107307H1-SSIT
16703-79	5102607H2-E	5107307H2-FF
16TPB002-05	5102607H2-EIVG	5107307H2-OO
4-FLEX	5102707H1	5107307H2-OOIT
5101507H9	5102707H2	5107307H2-SO
5101507H9--B	5102707H3	5107307H2-SOIT
5102107H1	5102807H1	5107307H2-SS
5102107H2	5102807H3	5107307H2-SSIT
5102107H2-E	5105307H9-FF-A	5107307H3-FF
5102507H1	5105307H9-OO-1	5107307H3-OO
5102507H1--IVG	5105307H9-OO-A	5107307H3-OOIT
5102507H1-E	5105307H9-SO-1	5107307H3-SO
5102507H1-ELE	5105307H9-SO-A	5107307H3-SOIT
5102507H1-X	5105307H9-SS-A	5107307H3-SS
5102507H2	5105307H9-SSVPA	5107307H3-SSIT
5102507H2--IVG	5105307H9-EOO-A	5152507T0--EA
5102507H2-E	5105307H9-ESO-A	5152507T1--C
5102507H2-EIVG	5105307H9-ESS-A	5152507T1--EA
5102507H2-ELE	5107307H1-FF	5152507T1-E
5102507H3	5107307H1-OO	5152507T1-E-C
5102507H3--IVG	5107307H1-OOIT	5152507T2--C
5102507H3-E	5107307H1-SO	5152507T2--EA
5102507H3-X	5107307H1-SOIT	5152507T2-E

5152507T2-E-C	5181507T7--B	5182107T1--A-K
5152507T3--EA	5181507T7--GPN	5182107T1--GA
5152507T3-E-C	5181507T7-E-B	5182107T1--GF
5152707T1-E	5181507T7-EGPN	5182107T1--LG
5152707T1-E-C	5181507T8--B	5182107T1-FULG
5152707T2-E	5181507T8--MDB	5182158T1--A-K
5152707T2-E-C	5181507T8-E-B	5182158T1-FULG
5152807T1-E	5181507T9--B	5182506T0--GA
5152807T1-E-C	5181507T9--EAB	5182506T1--GA
5152807T2-E	5181507T9-E-B	5182506T1-EGA
5152807T2-E-C	5181507T9-EK-B	5182506T2--GA
5172507T1--HR	5181507T9-V-B	5182506T3--GA
5172507T2--HR	5181507T9-25-B	51825070-TCA
5172507T3--HR	5181507T9-25BXB	5182507T0--GA
5181207T7--B	5181807T6	5182507T0--GAG
5181207T8--B	5181807T6--D	5182507T0-EGA
5181207T8--MTB	5181807T6--K	5182507T0-TCA
5181207T8--STB	5181807T6--EAA	5182507T0-TGA
5181207T9--B	5181807T7--B	5182507T0-TLE
5181207T9--MTB	5181807T7--KSH	5182507T0-VGA
5181307T4-OO-B	5181807T8--B	5182507T1--F2
5181307T4-SO-B	5181807T8-E-B	5182507T1--F3
5181307T4-SS-1	5181807T9--B	5182507T1--GA
5181307T4-SS-B	5182106T1-FULG	5182507T1--GAG
5181307T5-OO-D	5182107T0--A-K	5182507T1--LG
5181407T7--B	5182107T0--GF	5182507T1--LGG

5182507T1--SC	5182507T3--EGA	5182707T3--LG
5182507T1--EGA	5182507T3--ELG	5182707T3--FSLG
5182507T1--ELG	5182507T3--VGA	5182807T1--LG
5182507T1--ESG	5182558T1--GA	5182807T1--EKO
5182507T1--VGA	5182558T1--LG	5182807T1--ELG
5182507T1--XGA	5182558T3--LG	5182807T2--LG
5182507T1--XLG	5182603T2--LGE	5182807T2--EKO
5182507T1--FSLGG	5182606T3--GA	5182807T2--ELG
5182507T1A--LG	5182607T1--GA	5182807T3--LG
5182507T2--GA	5182607T2--LG	5185307T0--OO-1
5182507T2--GAG	5182607T2--LGE	5185307T0--SO-1
5182507T2--HR	5182607T2--EGA	5185307T0--SO-V
5182507T2--LG	5182607T3--GA	5185307T0--SS-V
5182507T2--SC	5182607T3--VGA	5185307T7--FF
5182507T2--EGA	5182658T1--LGE	5185307T7--FF-A
5182507T2--EGAG	5182658T2--LGE	5185307T7--FH
5182507T2--ELG	5182706T1--FULG	5185307T7--FH-A
5182507T2--VGA	5182707T0--GA	5185307T7--FO
5182507T2--XGA	5182707T1--GA	5185307T7--FO-A
5182507T2--FSLG	5182707T1--LG	5185307T7--HH-A
5182507T2B--GA	5182707T1--KFSG	5185307T7--NN-A
5182507T3--F2	5182707T1--FSLG	5185307T7--OO-1
5182507T3--GA	5182707T1--FULG	5185307T7--OO-A
5182507T3--GAG	5182707T2--GA	5185307T7--RR-A
5182507T3--LG	5182707T2--LG	5185307T7--SO-1
5182507T3--SC	5182707T2--FSLG	5185307T7--SO-A

5185307T7-SS-A	5185307T9-OOOAA	5195307A0G-SS-V
5185307T7-SSP1A	5185307T9-SO-1	5195307A7G-FF-1
5185307T7-WW-A	5185307T9-SO-A	5195307A7G-FF-A
5185307T8-FF-A	5185307T9-SS	5195307A7G-FO-1
5185307T8-NN-A	5185307T9-SS-A	5195307A7G-FO-A
5185307T8-OO	5185507T1--GA	5195307A7G-OO-1
5185307T8-OO-1	5187307T1-BB	5195307A7G-OO-A
5185307T8-OO-A	5187307T1-FF	5195307A7G-SO-1
5185307T8-OO-K	5187307T1-SB	5195307A7G-SO-A
5185307T8-OOOAA	5187307T1-SS	5195307A7G-SS-1
5185307T8-RR-A	5187307T2-SS	5195307A7G-SS-A
5185307T8-SO-1	5191307A4G-OO-1	5195307A8G-FF-1
5185307T8-SO-A	5191307A4G-OO-B	5195307A8G-FF-A
5185307T8-SS	5191307A4G-SO-1	5195307A8G-FO-1
5185307T8-SS-1	5191307A4G-SO-B	5195307A8G-FO-A
5185307T8-SS-A	5191307A4G-SS-1	5195307A8G-OO-1
5185307T8-SSCCA	5191307A4G-SS-B	5195307A8G-OO-A
5185307T8-WW-A	5195307A0G-FF-1	5195307A8G-SO-1
5185307T8-ZZ-A	5195307A0G-FF-V	5195307A8G-SO-A
5185307T9-FF-1	5195307A0G-FO-1	5195307A8G-SS-1
5185307T9-FF-A	5195307A0G-FO-V	5195307A8G-SS-A
5185307T9-FH-A	5195307A0G-OO-1	5195307A9G-FF-1
5185307T9-HF-A	5195307A0G-OO-V	5195307A9G-FF-A
5185307T9-NN-A	5195307A0G-SO-1	5195307A9G-FO-1
5185307T9-OO-1	5195307A0G-SO-V	5195307A9G-FO-A
5185307T9-OO-A	5195307A0G-SS-1	5195307A9G-OO-1

5195307A9G-OO-A	5202506A3	5202805A3
5195307A9G-SO-1	5202506A4	5202806A2
5195307A9G-SO-A	5202506F1--BE	5202806A3
5195307A9G-SS-1	5202506F2--BE	5202807D3
5195307A9G-SS-A	5202506F3--BE	5202807D3---G
5197307A1G-OO	5202506F3--BEC	5202B06F3--BE
5197307A1G-OOIT	5202506F4--BE	5202B06F3--BEC
5197307A1G-SO	5202507D3	5212107A3
5197307A1G-SOIT	5202507D4	5212507A1
5197307A1G-SS	5202566D3--BEC	5212507A2
5197307A1G-SSIT	5202605F2--BE	5212507A3
5197307A2G-OO	5202605F3--BE	5212507A3---G
5197307A2G-OOIT	5202606F1--BE	5212707A3
5197307A2G-SO	5202606F2--BE	5212807A3
5197307A2G-SOIT	5202606F3--BE	5212807A3--G
5197307A2G-SS	5202606F3--BEC	5212807A3-E
5197307A2G-SSIT	5202705A2	5231306D4-SS-B
5202106F1--BE	5202705A3	5231306D5-OO-B
5202505A1	5202705F1--BE	5231307H4-OO
5202505A2	5202705F2--BE	5231307H4-OO-B
5202505A3	5202705F3--BE	5231307H4-OOAEB
5202505A4	5202705F3--BEC	5231307H4-ODDPB
5202505F3--BE	5202706F1--BE	5231307H4-OOPTB
5202505F3--BEC	5202706F2--BE	5231307H4-SO-B
5202506A1	5202706F3--BE	5231307H4-SOAEB
5202506A2	5202706F3--BEC	5231307H4-SS-B

5231307H4-SSVPB	5231807H7--B	5232507H1--PH
5231307H5-NO-B	5231807H8--B	5232507H1-EIVG
5231307H5-NO-V	5231807H9--B	5232507H1-EP
5231307H5-OO	5232106D1--P	5232507H2--BH
5231307H5-OO-B	5232107H1--IVG	5232507H2--HB
5231307H5-OO-D	5232107H1--P	5232507H2--IVG
5231307H5-OO-V	5232107H2--P	5232507H2--P
5231307H5-ODPB	5232158H1--P	5232507H2-EIVG
5231307H5-OOPTB	5232506D0--P	5232507H2-EP
5231358G0-SS-V	5232506D1--IVG	5232507H3--IVG
5231358H0-SO-1	5232506D1--P	5232507H3--P
5231358H5-OO-D	5232506D1--P-H	5232507H3--PH
5231506D7--B	5232506D2--IVG	5232507H3-EIVG
5231506D8--B	5232506D2--P	5232507H3-EP
5231506D9	5232506D3--IVG	5232558H0--P
5231506D9--B	5232506D3--P	5232558H1--P
5231507H7	5232506D4--IVG	5232558H2--P
5231507H7--B	5232506D4--P	5232558H3--P
5231507H7-MRUM	5232507G1-PH-2O	5232606D2--P
5231507H8	5232507G2-PH-2O	5232607G1-PH-2O
5231507H8--B	5232507G3-PH-2O	5232607H1--P
5231507H8-V-B	5232507H0--P	5232607H2--P
5231507H9	5232507H0-VP	5232607H3--P
5231507H9--B	5232507H0-V-P	5232706D1--P
5231558H8--B	5232507H1--IVG	5232706D2--P
5231558H9--B	5232507H1--P	5232706D3--P

5232707G1--PH	5235306D0-SS-V	5235307H0-SO-1
5232707G2-PH-2O	5235306D7-FF-A	5235307H0-SO-V
5232707H1--P	5235306D7-OO-1	5235307H0-SS-V
5232707H2--P	5235306D7-OO-A	5235307H4-ODPB
5232707H3--P	5235306D7-ODPA	5235307H5-ODPB
5232758H1--P	5235306D7-ODPB	5235307H7-FF-A
5232758H2--P	5235306D7-SO-1	5235307H7-FH-A
5232758H3-EP	5235306D7-SO-A	5235307H7-FO-A
5232806D1--IVG	5235306D7-SS-A	5235307H7-HH-A
5232806D1--P	5235306D7-WW-A	5235307H7-NN-A
5232806D2--P	5235306D8-FF-A	5235307H7-OO
5232806D3--P	5235306D8-OO-1	5235307H7-OO-1
5232807H1--IVG	5235306D8-OO-A	5235307H7-OO-A
5232807H1--P	5235306D8-SO-1	5235307H7-ODPA
5232807H2--P	5235306D8-SO-A	5235307H7-SO
5232807H3--IVG	5235306D8-SS-A	5235307H7-SO-1
5232807H3--P	5235306D9-NN-A	5235307H7-SO-A
5232858H1--P	5235306D9-OO-1	5235307H7-SO-K
5232858H2--P	5235306D9-OO-A	5235307H7-SODPA
5232858H3--P	5235306D9-SO-1	5235307H7-SS-A
5232858H4--P	5235306D9-SO-A	5235307H7-SSDPA
5235306D0-FF-V	5235306D9-SS-A	5235307H7-WW-A
5235306D0-OO-1	5235307H0-FF-V	5235307H8-FF-A
5235306D0-OO-V	5235307H0-HH-V	5235307H8-NO-A
5235306D0-SO-1	5235307H0-OO-1	5235307H8-OO
5235306D0-SO-V	5235307H0-OO-V	5235307H8-OO-1

5235307H8-OO-A	5235358H0-OO-1	5237306D2-ODX
5235307H8-RO-A	5235358H0-SO-1	5237306D2-OOIT
5235307H8-RR-A	5235358H0-SO-V	5237306D2-SB
5235307H8-SO	5235358H0-SS-V	5237306D2-SO
5235307H8-SO-1	5235358H7-OO-1	5237306D2-SS
5235307H8-SO-A	5235358H7-OO-A	5237306D3-BB
5235307H8-SO-K	5235358H7-SO-1	5237306D3-FF
5235307H8-SODPA	5235358H7-SO-A	5237306D3-OO
5235307H8-SS	5235358H7-SS-A	5237306D3-SB
5235307H8-SS-A	5235358H8-OO-1	5237306D3-SO
5235307H8-SSDPA	5235358H8-OO-A	5237306D3-SS
5235307H8-WW-A	5235358H8-SO-1	5237307G3H-OO2O
5235307H8-ZO-A	5235358H8-SS-A	5237307G3H-SS2O
5235307H8-ZZ-A	5235358H9-OO-1	5237307H1-BB
5235307H9-FF	5235358H9-OO-A	5237307H1-FF
5235307H9-FF-A	5235358H9-SO-1	5237307H1-OO
5235307H9-NN-A	5235358H9-SS-A	5237307H1-OOIT
5235307H9-OO-1	5237306D1-BB	5237307H1-SB
5235307H9-OO-A	5237306D1-FF	5237307H1-SO
5235307H9-OOCEA	5237306D1-OO	5237307H1-SOIT
5235307H9-RR-A	5237306D1-SB	5237307H1-SS
5235307H9-SO-1	5237306D1-SO	5237307H1-SSDP
5235307H9-SO-A	5237306D1-SS	5237307H1-SSIT
5235307H9-SODPA	5237306D2-BB	5237307H2-BB
5235307H9-SS-A	5237306D2-FF	5237307H2-FF
5235307H9-SSDPA	5237306D2-OO	5237307H2-OO

5237307H2-OOIT	5237358H3-SO	5302505A3--SC
5237307H2-SB	5237358H3-SS	5302505A4
5237307H2-SO	5238306D1-OO	5302505A4--EA
5237307H2-SOIT	5238306D2-OO	5302505A4--KI
5237307H2-SS	5238306D3-OO	5302506A1
5237307H2-SSDP	5238307H1-OO	5302506A2
5237307H2-SSIT	5238307H1-SO	5302506A2--G
5237307H3-BB	5238307H1-SS	5302506A3
5237307H3-FF	5238307H1-SY	5302506A3--G
5237307H3-OO	5238307H1-YY	5302506A3--EA
5237307H3-ODDX	5238307H2-OO	5302506A3--KI
5237307H3-OOIT	5238307H2-SO	5302506A3--SC
5237307H3-SB	5238307H2-SS	5302506A4
5237307H3-SO	5238307H2-SY	5302506A4--G
5237307H3-SOIT	5238307H2-YY	5302506A4--EA
5237307H3-SS	5238307H3-OO	5302506A4--KI
5237307H3-SSIT	5238307H3-SO	5302706A3
5237358H1-FF	5238307H3-SS	5302805A3
5237358H1-OO	5238307H3-SY	5302805A3--G
5237358H1-SO	5238307H3-YY	5302805A3--EA
5237358H1-SS	5302105A1--A-K	5302806A1
5237358H2-FF	5302505A1	5302806A2
5237358H2-OO	5302505A2	5302806A3
5237358H2-SO	5302505A3	5302806A3--G
5237358H2-SS	5302505A3--EA	5302806A3--EA
5237358H3-OO	5302505A3--KI	5302806A4

5302806A4--EA	5401307A5-OO-B	5412507A4--EB
5312505A3	5402807A1	5412507A4--SC
5312505A3-E	5412107A1	5412707A1--EB
5312505A4	5412107A1--A-K	5412707A2
5312524A1--EB	5412107A1--FW	5412707A3
5312524A2	5412107A1--GF	5412807A1
5312524A2--EB	5412107A2	5412807A2
5312524A3	5412507A1	5412807A2--EA
5312524A3--G	5412507A1--EB	5412807A2--EB
5312524A3--EB	5412507A1--XEB	5412807A3
5312524A4	5412507A1-E	5412807A3--G
5312524A4--EB	5412507A1-XEB	5412807A3--EA
5312724A3	5412507A2	5412807A3--EB
5312805A3	5412507A2--EA	5412807A3-E
5312824A2	5412507A2--EB	5412807A3-EK
5312824A3	5412507A2-E	5412807A4
5312824A4	5412507A3	5415307A7--BRK
5382507H2	5412507A3--G	5422105A1--A-K
5382507H2--G	5412507A3--EA	5422105A1--GF
5382507H3	5412507A3--EB	5422106A1--A-K
5382507H3--G	5412507A3--SC	5422106A1--GF
5382507H4	5412507A3-E	5422505A1
5382507H4--G	5412507A3-V	5422505A1--EB
5382558K3	5412507A4	5422505A2
5382807H3	5412507A4--G	5422505A2--EA
5382858K3	5412507A4--EA	5422505A3

5422505A3--G	5422705A3--EB	5441307G4-SS-B
5422505A3--EA	5422706A3	5441307G4-SSVPB
5422505A3--EB	5422706A3--EB	5441307G4G-OO-1
5422505A3--FW	5422805A1	5441307G4G-OO-B
5422505A3--JK	5422805A2	5441307G4G-SO-1
5422505A3--KI	5422805A3	5441307G4G-SO-B
5422505A3--SC	5422805A3--G	5441307G4G-SS-1
5422505A3-E	5422805A3--EA	5441307G4G-SS-B
5422505A4	5422806A1	5441307G4G-SSVPB
5422505A4--EB	5422806A2	5441307H4-NO-B
5422505A4--KI	5422806A2--EA	5441307H4-NOIN
5422506A1	5422806A3	5441307H4-OO
5422506A1--EB	5422806A3--G	5441307H4-OO-B
5422506A2	5422806A3--EA	5441307H4-SO-B
5422506A2--EA	5422806A4	5441307H4-SS-B
5422506A2--EB	5441306A5-OOEIK	5441307H4-SSVP
5422506A2-E	5441306G4-OO-B	5441307H4-SSVPB
5422506A3	5441306G5-OO	5441307H4G-OO-1
5422506A3--G	5441306G5-OO-1	5441307H4G-OO-B
5422506A3--EA	5441306G5-OO-B	5441307H4G-OOOXQ
5422506A3--EB	5441307A4G-OO-B	5441307H4G-OOSD
5422506A3--SC	5441307A4G-SO-B	5441307H4G-OOTCB
5422506A4	5441307A4G-SS-B	5441307H4G-SO-1
5422506A4--EB	5441307G4-NO-B	5441307H4G-SO-B
5422506A4--SC	5441307G4-OO-B	5441307H4G-SOTC1
5422705A3	5441307G4-SO-B	5441307H4G-SOTCB

5441307H4G-SS-1	5441307I4G-SSVPB	5441507H7--B
5441307H4G-SS-B	5441307IA-OO-B	5441507H7--IOB
5441307H4G-SSVPB	5441358K4-OO-B	5441507H7-E-B
5441307H5-OO	5441358K4-SO-B	5441507H8
5441307H5-OO-1	5441358K4-SS-B	5441507H8--B
5441307H5-OO-B	5441358K4G-OO-1	5441507H9
5441307H5-OOPIB	5441358K4G-OO-B	5441507H9--B
5441307H5G-OO-1	5441358K4G-SO-1	5441507H9-E-B
5441307H5G-OO-B	5441358K4G-SO-B	5441507H9-25-B
5441307H5G-OOSD	5441358K5-OO	5441558K7--B
5441307H5G-OOSD1	5441358K5-OO-B	5441558K8--B
5441307H5G-OOSDK	5441358K8G-SS-B	5441558K9
5441307H5G-OOTCB	5441358K9G-SS-A	5441558K9--B
5441307H9G-OO-A	5441358M4-OO-B	5441558M8-25-B
5441307I4-NO-B	5441358M4-SO-B	5441558M9-25-B
5441307I4-OO-B	5441358M4-SS-B	5441807H7-E-B
5441307I4-SO-B	5441358M4G-OO-1	5441807H8--B
5441307I4-SS-B	5441358M4G-OO-B	5441807H9--B
5441307I4-SSVP	5441358M4G-SO-1	5442106G1
5441307I4-SSVPB	5441358M4G-SO-B	5442107H1
5441307I4G-OO-1	5441358M4G-SS-1	5442107H2
5441307I4G-OO-B	5441358M4G-SS-B	5442107H3
5441307I4G-SO-1	5441506G7--B	5442158K1--GF
5441307I4G-SO-B	5441506G8--B	5442506G0
5441307I4G-SS-1	5441506G9--B	5442506G1
5441307I4G-SS-B	5441507H7	5442506G1--IVG

5442506G2	5442507H1--HB	5442558K2
5442506G2--IVG	5442507H1--IVG	5442558K2-E
5442506G2-E	5442507H1--SI	5442558K3
5442506G3	5442507H1-E	5442558K3--IVG
5442506G3--IVG	5442507H2	5442558K3-E
5442506G4	5442507H2--G	5442558M0
5442507A1--HF	5442507H2--HB	5442558M1
5442507A1--HFG	5442507H2--IVG	5442558M1--EB
5442507A2--HF	5442507H2-E	5442558M2
5442507A2--HFG	5442507H3	5442558M2--EB
5442507A3--HF	5442507H3--G	5442558M3
5442507A3--HFG	5442507H3--AO	5442558M3--EB
5442507G0	5442507H3--IVG	5442558M4
5442507G1	5442507H3-E	5442606G2-EAB
5442507G1--G	5442507H3-EIVG	5442607G2--ABG
5442507G2	5442507H4	5442607H1
5442507G2--G	5442507I1	5442706G1
5442507G2-E	5442507I1-E	5442706G2
5442507G3	5442507I2	5442706G3
5442507G3--G	5442507I2-E	5442706G4
5442507H0	5442507I3	5442707H1
5442507H0--G	5442507I3-E	5442707H1-VSI
5442507H0--HB	5442558K0	5442707H2
5442507H0-E	5442558K1	5442707H2--IVG
5442507H1	5442558K1--IVG	5442707H3
5442507H1--G	5442558K1-E	5442758K1

5442758K2	5445306G7-SS-A	5445306G9G-SO-1
5442806G3--IVG	5445306G7G-OO-1	5445306G9G-SO-A
5442807H1	5445306G7G-OO-A	5445306G9G-SS-1
5442807H1--IVG	5445306G7G-SO-1	5445306G9G-SS-A
5442807H2	5445306G7G-SO-A	5445307A0G-OO-V
5442807H3	5445306G7G-SS-1	5445307A0G-SO-V
5442858K1	5445306G7G-SS-A	5445307A0G-SS-V
5442858K2	5445306G8-FF-A	5445307A7G-OO-A
5442858K3	5445306G8-HH-A	5445307A7G-SO-A
5445306A7-HHEIK	5445306G8-OO-A	5445307A7G-SS-A
5445306G0-FF-V	5445306G8-SO-A	5445307A8G-OO-A
5445306G0-OO-V	5445306G8-SS-1	5445307A8G-SO-A
5445306G0-SO-V	5445306G8-SS-A	5445307A8G-SS-A
5445306G0-SS-V	5445306G8G-OO-1	5445307A9G-OO-A
5445306G0G-OO-1	5445306G8G-OO-A	5445307A9G-SO-A
5445306G0G-OO-V	5445306G8G-SO-1	5445307A9G-SS-A
5445306G0G-SO-1	5445306G8G-SO-A	5445307G0-FF-V
5445306G0G-SO-V	5445306G8G-SS-1	5445307G0-FO-V
5445306G0G-SS-1	5445306G8G-SS-A	5445307G0-OO-V
5445306G0G-SS-V	5445306G9-FF-A	5445307G0-SO-V
5445306G7-FF-A	5445306G9-OO-A	5445307G0-SS-V
5445306G7-FH-A	5445306G9-SO	5445307G0G-OO-1
5445306G7-FO-A	5445306G9-SO-A	5445307G0G-OO-V
5445306G7-HH-A	5445306G9-SS-A	5445307G0G-SO-1
5445306G7-OO-A	5445306G9G-OO-1	5445307G0G-SO-V
5445306G7-SO-A	5445306G9G-OO-A	5445307G0G-SS-1

5445307G0G-SS-V	5445307G9-SO-A	5445307H7-SO-A
5445307G7-FF-A	5445307G9-SS-A	5445307H7-SS-A
5445307G7-FO-A	5445307G9G-OO-1	5445307H7-EOO-A
5445307G7-OO-A	5445307G9G-OO-A	5445307H7-ESS-A
5445307G7-SO-A	5445307G9G-SO-1	5445307H7G-FF-1
5445307G7-SS-A	5445307G9G-SO-A	5445307H7G-FF-A
5445307G7G-OO-1	5445307G9G-SS-1	5445307H7G-FH-A
5445307G7G-OO-A	5445307G9G-SS-A	5445307H7G-FO-A
5445307G7G-SO-1	5445307H0-FF-V	5445307H7G-HH-1
5445307G7G-SO-A	5445307H0-OO-V	5445307H7G-HH-A
5445307G7G-SS-1	5445307H0-SF-V	5445307H7G-OO-1
5445307G7G-SS-A	5445307H0-SO-V	5445307H7G-OO-A
5445307G8-FF-A	5445307H0-SS-V	5445307H7G-SO-1
5445307G8-FO-A	5445307H0G-OO	5445307H7G-SO-A
5445307G8-OO-A	5445307H0G-OO-1	5445307H7G-SS-1
5445307G8-SO-A	5445307H0G-OO-V	5445307H7G-SS-A
5445307G8-SS-A	5445307H0G-SO-V	5445307H8-FF-A
5445307G8G-OO-1	5445307H0G-SO-1	5445307H8-OO-A
5445307G8G-OO-A	5445307H0G-SO-V	5445307H8-SF-A
5445307G8G-SO-1	5445307H0G-SS-1	5445307H8-SO-A
5445307G8G-SO-A	5445307H0G-SS-V	5445307H8-SS-A
5445307G8G-SS-1	5445307H7-FF-A	5445307H8-WW-A
5445307G8G-SS-A	5445307H7-FH-A	5445307H8-EOO-A
5445307G9-FF-A	5445307H7-FO-A	5445307H8-ESO-A
5445307G9-FO-A	5445307H7-HH-A	5445307H8-ESS-A
5445307G9-OO-A	5445307H7-OO-A	5445307H8G-OO-1

5445307H8G-OO-A	5445307I0G-OO-1	5445307I8G-FF-1
5445307H8G-SO-1	5445307I0G-OO-V	5445307I8G-FF-A
5445307H8G-SO-A	5445307I0G-SO-1	5445307I8G-FO-1
5445307H8G-SS-1	5445307I0G-SO-V	5445307I8G-FO-A
5445307H8G-SS-A	5445307I0G-SS-1	5445307I8G-OO-1
5445307H9-FF-A	5445307I0G-SS-V	5445307I8G-OO-A
5445307H9-OO-A	5445307I7-FF-A	5445307I8G-SO-1
5445307H9-SO-A	5445307I7-FO-A	5445307I8G-SO-A
5445307H9-SS-A	5445307I7-OO-A	5445307I8G-SS-1
5445307H9G-OO-1	5445307I7-SO-A	5445307I8G-SS-A
5445307H9G-OO-A	5445307I7-SS-A	5445307I9-FF-A
5445307H9G-SO-1	5445307I7G-FF-1	5445307I9-FO-A
5445307H9G-SO-A	5445307I7G-FF-A	5445307I9-OO-A
5445307H9G-SS-1	5445307I7G-FO-1	5445307I9-SO-A
5445307H9G-SS-A	5445307I7G-FO-A	5445307I9-SS-A
5445307H0G-SO	5445307I7G-OO-1	5445307I9G-FF-1
5445307H0G-SS	5445307I7G-OO-A	5445307I9G-FF-A
5445307I0-FF-V	5445307I7G-SO-1	5445307I9G-FO-1
5445307I0-FO-V	5445307I7G-SO-A	5445307I9G-FO-A
5445307I0-OO-V	5445307I7G-SS-1	5445307I9G-OO-1
5445307I0-SO-V	5445307I7G-SS-A	5445307I9G-OO-A
5445307I0-SS-V	5445307I8-FF-A	5445307I9G-SO-1
5445307I0G-FF-1	5445307I8-FO-A	5445307I9G-SO-A
5445307I0G-FF-V	5445307I8-OO-A	5445307I9G-SS-1
5445307I0G-FO-1	5445307I8-SO-A	5445307I9G-SS-A
5445307I0G-FO-V	5445307I8-SS-A	5445358K0-FF-V

5445358K0-OO-V	5445358K7G-SO-1	5445358KOG-OO
5445358K0-SO-V	5445358K7G-SO-A	5445358M0-FF-V
5445358K0-SS-V	5445358K7G-SS-1	5445358M0-FO-V
5445358KOG-OO-1	5445358K7G-SS-A	5445358M0-OO-V
5445358KOG-OO-V	5445358K8-FF-A	5445358M0-SO-V
5445358KOG-SO	5445358K8-OO-A	5445358M0-SS-V
5445358KOG-SO-1	5445358K8-SO-A	5445358M0-OO-V
5445358KOG-SO-V	5445358K8-SS-A	5445358M0-SO-V
5445358KOG-SS	5445358K8G-OO-1	5445358M0-SS-V
5445358KOG-SS-1	5445358K8G-OO-A	5445358M0G-FF-1
5445358KOG-SS-V	5445358K8G-SO-1	5445358M0G-FF-V
5445358K7-FF-A	5445358K8G-SO-A	5445358M0G-FO-1
5445358K7-FH-A	5445358K8G-SS-1	5445358M0G-FO-V
5445358K7-FO-A	5445358K8G-SS-A	5445358M0G-OO-1
5445358K7-HH-1	5445358K9-FF	5445358M0G-OO-V
5445358K7-HH-A	5445358K9-FF-A	5445358M0G-SO-1
5445358K7-OO-A	5445358K9-OO-1	5445358M0G-SO-V
5445358K7-SO-A	5445358K9-OO-A	5445358M0G-SS-1
5445358K7-SS-A	5445358K9-SO-A	5445358M0G-SS-V
5445358K7G-FF-1	5445358K9-SS-A	5445358M7-FF-A
5445358K7G-FF-A	5445358K9G-OO-1	5445358M7-FO-A
5445358K7G-FH-A	5445358K9G-OO-A	5445358M7-OO-A
5445358K7G-FO-A	5445358K9G-SO-1	5445358M7-SO-A
5445358K7G-HH-A	5445358K9G-SO-A	5445358M7-SS-A
5445358K7G-OO-1	5445358K9G-SS-1	5445358M7-FF-A
5445358K7G-OO-A	5445358K9G-SS-A	5445358M7-OO-A

5445358M7-SO-A	5445358M8G-SO-A	5447306G1G-OO
5445358M7-SS-A	5445358M8G-SS-1	5447306G1G-SO
5445358M7G-FF-1	5445358M8G-SS-A	5447306G1G-SS
5445358M7G-FF-A	5445358M9-FF-A	5447306G2-FF
5445358M7G-FO-1	5445358M9-FO-A	5447306G2-OO
5445358M7G-OO-1	5445358M9-OO-A	5447306G2-SO
5445358M7G-OO-A	5445358M9-SO-A	5447306G2-SS
5445358M7G-SO-1	5445358M9-SS-A	5447306G2G-OO
5445358M7G-SO-A	5445358M9-OO-A	5447306G2G-SO
5445358M7G-SS-1	5445358M9-SO-A	5447306G2G-SS
5445358M7G-SS-A	5445358M9-SS-A	5447306G3-FF
5445358M8-FF-A	5445358M9G-FF-1	5447306G3-OO
5445358M8-FO-A	5445358M9G-FF-A	5447306G3-SO
5445358M8-OO-A	5445358M9G-FO-1	5447306G3-SS
5445358M8-SO-A	5445358M9G-FO-A	5447306G3-SSAC
5445358M8-SS-A	5445358M9G-OO-1	5447306G3G-OO
5445358M8-OO-A	5445358M9G-OO-A	5447306G3G-SO
5445358M8-SO-A	5445358M9G-SO-1	5447306G3G-SS
5445358M8-SS-A	5445358M9G-SO-A	5447307G1-FF
5445358M8G-FF-1	5445358M9G-SS-1	5447307G1-OO
5445358M8G-FF-A	5445358M9G-SS-A	5447307G1-OOIT
5445358M8G-FO-1	5447306G1-FF	5447307G1-SO
5445358M8G-FO-A	5447306G1-OO	5447307G1-SOIT
5445358M8G-OO-1	5447306G1-SO	5447307G1-SS
5445358M8G-OO-A	5447306G1-SS	5447307G1-SSAC
5445358M8G-SO-1	5447306G1G-SSAC	5447307G1-SSIT

5447307G1G-OO	5447307G3-SSIT	5447307H2-OOIT
5447307G1G-OOIT	5447307G3G-OO	5447307H2-SO
5447307G1G-SO	5447307G3G-OOIT	5447307H2-SOHB
5447307G1G-SOIT	5447307G3G-SO	5447307H2-SOIT
5447307G1G-SS	5447307G3G-SOIT	5447307H2-SS
5447307G1G-SSIT	5447307G3G-SS	5447307H2-SSAC
5447307G2-FF	5447307G3G-SSIT	5447307H2-SSHB
5447307G2-OO	5447307H1-FF	5447307H2-SSIT
5447307G2-OOIT	5447307H1-OO	5447307H2G-OO
5447307G2-SO	5447307H1-OOAC	5447307H2G-SO
5447307G2-SOIT	5447307H1-OOIT	5447307H2G-SOIT
5447307G2-SS	5447307H1-SF	5447307H2G-SS
5447307G2-SSAC	5447307H1-SO	5447307H2G-SSIT
5447307G2-SSIT	5447307H1-SOIT	5447307H3-FF
5447307G2G-OO	5447307H1-SS	5447307H3-OO
5447307G2G-OOIT	5447307H1-SSAC	5447307H3-OOAC
5447307G2G-SO	5447307H1-SSIT	5447307H3-OOIT
5447307G2G-SOIT	5447307H1G-FO	5447307H3-SO
5447307G2G-SS	5447307H1G-OO	5447307H3-SOIT
5447307G2G-SSIT	5447307H1G-SO	5447307H3-SS
5447307G3-OO	5447307H1G-SOIT	5447307H3-SSAC
5447307G3-OOIT	5447307H1G-SS	5447307H3-SSIT
5447307G3-SO	5447307H1G-SSIT	5447307H3G-OO
5447307G3-SOIT	5447307H2-FF	5447307H3G-SO
5447307G3-SS	5447307H2-OO	5447307H3G-SOIT
5447307G3-SSAC	5447307H2-OOAC	5447307H3G-SS

5447307H3G-SSIT	5447307I2G-SO	5447358K1G-FO
5447307H3GMSS	5447307I2G-SOIT	5447358K1G-OO
5447307I1-OO	5447307I2G-SS	5447358K1G-SO
5447307I1-OOAC	5447307I2G-SSIT	5447358K1G-SOIT
5447307I1-OOIT	5447307I3-OO	5447358K1G-SS
5447307I1-SO	5447307I3-OOAC	5447358K1G-SSIT
5447307I1-SOIT	5447307I3-OOIT	5447358K2-FF
5447307I1-SS	5447307I3-SO	5447358K2-OO
5447307I1-SSAC	5447307I3-SOIT	5447358K2-OOIT
5447307I1-SSIT	5447307I3-SS	5447358K2-SO
5447307I1G-OO	5447307I3-SSAC	5447358K2-SOIT
5447307I1G-OOIT	5447307I3-SSIT	5447358K2-SS
5447307I1G-SO	5447307I3G-OO	5447358K2-SSAC
5447307I1G-SOIT	5447307I3G-OOIT	5447358K2-SSIT
5447307I1G-SS	5447307I3G-SO	5447358K2G-FO
5447307I1G-SSIT	5447307I3G-SOIT	5447358K2G-SO
5447307I2-OO	5447307I3G-SS	5447358K2G-SOIT
5447307I2-OOAC	5447307I3G-SSIT	5447358K2G-SS
5447307I2-OOIT	5447358K1-FF	5447358K2G-SSIT
5447307I2-SO	5447358K1-OO	5447358K3-FF
5447307I2-SOIT	5447358K1-OOIT	5447358K3-OO
5447307I2-SS	5447358K1-SO	5447358K3-OOIT
5447307I2-SSAC	5447358K1-SOIT	5447358K3-SO
5447307I2-SSIT	5447358K1-SS	5447358K3-SOIT
5447307I2G-OO	5447358K1-SSAC	5447358K3-SS
5447307I2G-OOIT	5447358K1-SSIT	5447358K3-SSIT

5447358K3G-SO	5447358M3G-SO	5448307G1G-SY
5447358K3G-SS	5447358M3G-SOIT	5448307G1G-YY
5447358K3G-SSIT	5447358M3G-SS	5448307G2G-OO
5447358M1-OO	5447358M3G-SSIT	5448307G2G-SO
5447358M1-SO	5448306G1G-OO	5448307G2G-SS
5447358M1-SS	5448306G1G-SO	5448307G2G-SY
5447358M1G-OO	5448306G1G-SS	5448307G2G-YY
5447358M1G-OOIT	5448306G2G-OO	5448307G3G-OO
5447358M1G-SO	5448306G2G-SO	5448307G3G-SO
5447358M1G-SOIT	5448306G2G-SS	5448307G3G-SS
5447358M1G-SS	5448306G3G-OO	5448307G3G-SY
5447358M1G-SSIT	5448306G3G-SO	5448307G3G-YY
5447358M2-OO	5448306G3G-SS	5448307H1G-OO
5447358M2-SO	5448307A1G-OO	5448307H1G-SO
5447358M2-SS	5448307A1G-SO	5448307H1G-SS
5447358M2G-OO	5448307A1G-SS	5448307H1G-SSAC
5447358M2G-OOIT	5448307A2G-OO	5448307H1G-SSPZ
5447358M2G-SO	5448307A2G-SO	5448307H1G-SY
5447358M2G-SOIT	5448307A2G-SS	5448307H1G-YY
5447358M2G-SS	5448307A3G-OO	5448307H2G-OO
5447358M2G-SSIT	5448307A3G-SO	5448307H2G-SO
5447358M3-OO	5448307A3G-SS	5448307H2G-SS
5447358M3-SO	5448307G1G-OO	5448307H2G-SSPZ
5447358M3-SS	5448307G1G-OOAC	5448307H2G-SY
5447358M3G-OO	5448307G1G-SO	5448307H2G-YY
5447358M3G-OOIT	5448307G1G-SS	5448307H3G-OO

5448307H3G-SO	5448358K3G-SS	5462507G3
5448307H3G-SS	5448358K3G-SY	5462558M1
5448307H3G-SSPZ	5448358K3G-YY	5462558M2
5448307H3G-SY	5448358M1G-OO	5462558M3
5448307H3G-YY	5448358M1G-SO	5465307G0G-FF-1
5448307I1G-OO	5448358M1G-SS	5465307G0G-FF-B
5448307I1G-SO	5448358M1G-SY	5465307G0G-FF-V
5448307I1G-SS	5448358M1G-YY	5465307G0G-FO-1
5448307I2G-OO	5448358M2G-OO	5465307G0G-FO-B
5448307I2G-SO	5448358M2G-SO	5465307G0G-FO-V
5448307I2G-SS	5448358M2G-SS	5465307G0G-OO-1
5448307I3G-OO	5448358M2G-SY	5465307G0G-OO-B
5448307I3G-SO	5448358M2G-YY	5465307G0G-OO-V
5448307I3G-SS	5448358M3G-OO	5465307G0G-SO-1
5448358K1G-OO	5448358M3G-SO	5465307G0G-SO-B
5448358K1G-SO	5448358M3G-SS	5465307G0G-SO-V
5448358K1G-SS	5448358M3G-SY	5465307G0G-SS-1
5448358K1G-SY	5448358M3G-YY	5465307G0G-SS-B
5448358K1G-YY	5461307G4-OO-B	5465307G0G-SS-V
5448358K2G-OO	5461307G4-SO-B	5465307G7G-OO-A
5448358K2G-SO	5461307G4-SS-B	5465307G8G-FF-1
5448358K2G-SS	5461358M4-OO-B	5465307G8G-FF-A
5448358K2G-SY	5461358M4-SO-B	5465307G8G-FO-1
5448358K2G-YY	5461358M4-SS-B	5465307G8G-FO-A
5448358K3G-OO	5462507G1	5465307G8G-OO-1
5448358K3G-SO	5462507G2	5465307G8G-OO-A

5465307G8G-SO-1	5465358M0G-SS-1	5467307G1G-SO
5465307G8G-SO-A	5465358M0G-SS-B	5467307G1G-SOIT
5465307G8G-SS-1	5465358M0G-SS-V	5467307G1G-SS
5465307G8G-SS-A	5465358M7G-SO-A	5467307G1G-SSIT
5465307G9G-FF-1	5465358M8G-FF-1	5467307G2G-OO
5465307G9G-FF-A	5465358M8G-FF-A	5467307G2G-OOIT
5465307G9G-FO-1	5465358M8G-FO-1	5467307G2G-SO
5465307G9G-FO-A	5465358M8G-FO-A	5467307G2G-SOIT
5465307G9G-OO-1	5465358M8G-OO-1	5467307G2G-SS
5465307G9G-OO-A	5465358M8G-OO-A	5467307G2G-SSIT
5465307G9G-SO-1	5465358M8G-SO-1	5467307G3G-OO
5465307G9G-SO-A	5465358M8G-SO-A	5467307G3G-OOIT
5465307G9G-SS-1	5465358M8G-SS-1	5467307G3G-SO
5465307G9G-SS-A	5465358M8G-SS-A	5467307G3G-SOIT
5465358M0G-FF-1	5465358M9G-FF-1	5467307G3G-SS
5465358M0G-FF-B	5465358M9G-FF-A	5467307G3G-SSIT
5465358M0G-FF-V	5465358M9G-FO-1	5467358M1G-OO
5465358M0G-FO-1	5465358M9G-FO-A	5467358M1G-SO
5465358M0G-FO-B	5465358M9G-OO-1	5467358M1G-SS
5465358M0G-FO-V	5465358M9G-OO-A	5467358M2G-OO
5465358M0G-OO-1	5465358M9G-SO-1	5467358M2G-SO
5465358M0G-OO-B	5465358M9G-SO-A	5467358M2G-SS
5465358M0G-OO-V	5465358M9G-SS-1	5467358M3G-OO
5465358M0G-SO-1	5465358M9G-SS-A	5467358M3G-SO
5465358M0G-SO-B	5467307G1G-OO	5467358M3G-SS
5465358M0G-SO-V	5467307G1G-OOIT	5468307G1

5468307G1G-OO	5468358M3G-SO	5472507F4-E
5468307G1G-SO	5468358M3G-SS	5472558G1
5468307G1G-SS	5468358M3G-SY	5472558G2
5468307G1G-SY	5468358M3G-YY	5472558G3
5468307G1G-YY	5471307F4-OO-B	5475307F0-OO-V
5468307G2G-OO	5471307F4-SO-B	5475307F0-SO-V
5468307G2G-SO	5471307F4-SS-B	5475307F0-SS-V
5468307G2G-SS	5471307F4G-OO-B	5475307F0G-OO-1
5468307G2G-SY	5471307F4G-SS-1	5475307F0G-OO-V
5468307G2G-YY	5471307F4G-SS-B	5475307F0G-SO-1
5468307G3G-OO	5471358G4-OO-B	5475307F0G-SO-V
5468307G3G-SO	5471358G4-SO-B	5475307F0G-SS-1
5468307G3G-SS	5471358G4-SS	5475307F0G-SS-V
5468307G3G-SY	5471358G4-SS-B	5475307F7-OO-A
5468307G3G-YY	5471358G4G-OO-B	5475307F7-SO-A
5468358M1G-OO	5471358G4G-SS-1	5475307F7-SS-A
5468358M1G-SO	5471358G4G-SS-B	5475307F7G-OO-1
5468358M1G-SS	5471558G0-25-V	5475307F7G-OO-A
5468358M1G-SY	5471558G9-25-B	5475307F7G-SO-1
5468358M1G-YY	5472507F1	5475307F7G-SO-A
5468358M2G-OO	5472507F1--G	5475307F7G-SS-1
5468358M2G-SO	5472507F2	5475307F7G-SS-A
5468358M2G-SS	5472507F2--G	5475307F8-OO-A
5468358M2G-SY	5472507F2-E	5475307F8-SO-A
5468358M2G-YY	5472507F3	5475307F8-SS-A
5468358M3G-OO	5472507F3-E	5475307F8G-OO-1

5475307F8G-OO-A	5475358G7G-OO-1	5475358G9G-SS-A
5475307F8G-SO-1	5475358G7G-OO-A	5477307F1-OO
5475307F8G-SO-A	5475358G7G-SO-1	5477307F1-SO
5475307F8G-SS-1	5475358G7G-SO-A	5477307F1-SS
5475307F8G-SS-A	5475358G7G-SS-1	5477307F1-SSAC
5475307F9-OO-A	5475358G7G-SS-A	5477307F1G-OO
5475307F9-SO-A	5475358G8-FF-A	5477307F1G-OOIT
5475307F9-SS-A	5475358G8-OO-A	5477307F1G-SO
5475307F9G-OO-1	5475358G8-SO-A	5477307F1G-SOIT
5475307F9G-OO-A	5475358G8-SS-A	5477307F1G-SS
5475307F9G-SO-1	5475358G8G-OO-1	5477307F1G-SSIT
5475307F9G-SO-A	5475358G8G-OO-A	5477307F2-OO
5475307F9G-SS-1	5475358G8G-SO-1	5477307F2-SO
5475307F9G-SS-A	5475358G8G-SO-A	5477307F2-SS
5475358G0-OO-V	5475358G8G-SS-1	5477307F2-SSAC
5475358G0-SO-V	5475358G8G-SS-A	5477307F2G-OO
5475358G0-SS-V	5475358G9-OO-1	5477307F2G-OOIT
5475358G0G-OO-1	5475358G9-OO-A	5477307F2G-SO
5475358G0G-OO-V	5475358G9-SO-A	5477307F2G-SOIT
5475358G0G-SO-1	5475358G9-SS-1	5477307F2G-SS
5475358G0G-SO-V	5475358G9-SS-A	5477307F2G-SSIT
5475358G0G-SS-1	5475358G9G-OO-1	5477307F3-OO
5475358G0G-SS-V	5475358G9G-OO-A	5477307F3-SO
5475358G7-OO-A	5475358G9G-SO-1	5477307F3-SS
5475358G7-SO-A	5475358G9G-SO-A	5477307F3-SSAC
5475358G7-SS-A	5475358G9G-SS-1	5477307F3G-OO

5477307F3G-OOIT	5477358G3G-OO	5478358G2G-OO
5477307F3G-SO	5477358G3G-OOIT	5478358G2G-SO
5477307F3G-SOIT	5477358G3G-SO	5478358G2G-SS
5477307F3G-SS	5477358G3G-SOIT	5478358G2G-SY
5477307F3G-SSIT	5477358G3G-SS	5478358G2G-YY
5477358G1-OO	5477358G3G-SSIT	5478358G3G-OO
5477358G1-SO	5478307F1G-OO	5478358G3G-SO
5477358G1-SS	5478307F1G-SO	5478358G3G-SS
5477358G1G-OO	5478307F1G-SS	5478358G3G-SY
5477358G1G-OOIT	5478307F1G-SY	5478358G3G-YY
5477358G1G-SO	5478307F1G-YY	5481307G4-SS-B
5477358G1G-SOIT	5478307F2G-OO	5481358M4-SS-B
5477358G1G-SS	5478307F2G-SO	5482507G1
5477358G1G-SSIT	5478307F2G-SS	5482507G2
5477358G2-OO	5478307F2G-SY	5482507G3
5477358G2-SO	5478307F2G-YY	5482507G3--G
5477358G2-SS	5478307F3G-OO	5482507G4
5477358G2G-OO	5478307F3G-SO	5482558M1
5477358G2G-OOIT	5478307F3G-SS	5482558M2
5477358G2G-SO	5478307F3G-SY	5482558M3
5477358G2G-SOIT	5478307F3G-YY	5485307G0-FF-V
5477358G2G-SS	5478358G1G-OO	5485307G0-FO-V
5477358G2G-SSIT	5478358G1G-SO	5485307G0-OO-1
5477358G3-OO	5478358G1G-SS	5485307G0-OO-V
5477358G3-SO	5478358G1G-SY	5485307G0-SO-1
5477358G3-SS	5478358G1G-YY	5485307G0-SO-V

5485307G0-SS-V	5485358M8-OO-1	5487307G3-SS
5485307G7-OO-A	5485358M8-OO-A	5487358M1-BB
5485307G7-SO-A	5485358M8-SO-1	5487358M1-OO
5485307G8-FF-A	5485358M8-SO-A	5487358M1-SB
5485307G8-FO-A	5485358M8-SS-A	5487358M1-SO
5485307G8-OO-1	5485358M9-FF-A	5487358M1-SS
5485307G8-OO-A	5485358M9-FO-A	5487358M2-BB
5485307G8-SO-1	5485358M9-OO-1	5487358M2-OO
5485307G8-SO-A	5485358M9-OO-A	5487358M2-SB
5485307G8-SS-A	5485358M9-SO-1	5487358M2-SO
5485307G9-FF-A	5485358M9-SO-A	5487358M2-SS
5485307G9-FO-A	5485358M9-SS-A	5487358M3-BB
5485307G9-OO-1	5487307G1-BB	5487358M3-OO
5485307G9-OO-A	5487307G1-OO	5487358M3-SB
5485307G9-SO-1	5487307G1-SB	5487358M3-SO
5485307G9-SO-A	5487307G1-SO	5487358M3-SS
5485307G9-SS-A	5487307G1-SS	5488307G1-OO
5485358M0-FF-V	5487307G2-BB	5488307G1-SO
5485358M0-FO-V	5487307G2-OO	5488307G1-SS
5485358M0-OO-1	5487307G2-SB	5488307G1-SY
5485358M0-OO-V	5487307G2-SO	5488307G1-YY
5485358M0-SO-1	5487307G2-SS	5488307G2-OO
5485358M0-SO-V	5487307G3-BB	5488307G2-SO
5485358M0-SS-V	5487307G3-OO	5488307G2-SS
5485358M8-FF-A	5487307G3-SB	5488307G2-SY
5485358M8-FO-A	5487307G3-SO	5488307G2-YY

5488307G3-OO	5491307H4-SOHBB	5492507H1-S
5488307G3-SO	5491307H4-SS-B	5492507H2
5488307G3-SS	5491307H4-SSHB1	5492507H2--HB
5488307G3-SY	5491307H4-SSHBB	5492507H2-S
5488307G3-YY	5491307H4-SSVPB	5492507H3
5488358M1-OO	5491307H4G-NO-1	5492507H3--HB
5488358M1-SO	5491307H4G-NO-B	5492507H3-S
5488358M1-SS	5491307H4G-OO-1	549307H8-OO-A
5488358M1-SY	5491307H4G-OO-B	549307H9-OO-A
5488358M1-YY	5491307H4G-OOHB1	5495307H0-FF-V
5488358M2-OO	5491307H4G-OOHBB	5495307H0-OO-V
5488358M2-SO	5491307H4G-SO-1	5495307H0-OOHBV
5488358M2-SS	5491307H4G-SO-B	5495307H0-SO-V
5488358M2-SY	5491307H4G-SOHB1	5495307H0-SOHBV
5488358M2-YY	5491307H4G-SOHBB	5495307H0-SS-V
5488358M3-OO	5491307H4G-SS-1	5495307H0-SSHBV
5488358M3-SO	5491307H4G-SS-B	5495307H0G-OO-1
5488358M3-SS	5491307H4G-SSHB1	5495307H0G-OO-V
5488358M3-SY	5491307H4G-SSHBB	5495307H0G-OOHB1
5488358M3-YY	5491307H4G-SSVPB	5495307H0G-OOHBV
5491307H4-NO-B	5491307HC-FF-V	5495307H0G-SO-1
5491307H4-OO-B	5491307HC-OO-V	5495307H0G-SO-V
5491307H4-OOHB1	5491307HC-OO-V	5495307H0G-SOHB1
5491307H4-OOHBB	5491507H9-25-B	5495307H0G-SOHBV
5491307H4-SO-B	5492507H1	5495307H0G-SS-1
5491307H4-SOHB1	5492507H1--HB	5495307H0G-SS-V

5495307H0G-SSHB1	5495307H7G-SS-1	5495307H9-SO-A
5495307H0G-SSHBV	5495307H7G-SS-A	5495307H9-SOHBA
5495307H7-FF-A	5495307H7G-SSHB1	5495307H9-SS-A
5495307H7-HH-A	5495307H7G-SSHBA	5495307H9-SSHBA
5495307H7-OO-A	5495307H8-FF-A	5495307H9G-OO-1
5495307H7-OOHBA	5495307H8-OO-A	5495307H9G-OO-A
5495307H7-SO-A	5495307H8-OOHBA	5495307H9G-OOHB1
5495307H7-SOHBA	5495307H8-SO-A	5495307H9G-OOHBA
5495307H7-SS-A	5495307H8-SOHBA	5495307H9G-SO-1
5495307H7-SSHBA	5495307H8-SS-A	5495307H9G-SO-A
5495307H7G-FF-1	5495307H8-SSHBA	5495307H9G-SOHB1
5495307H7G-FF-A	5495307H8G-OO-1	5495307H9G-SOHBA
5495307H7G-FH-1	5495307H8G-OO-A	5495307H9G-SS-1
5495307H7G-FH-A	5495307H8G-OOHB1	5495307H9G-SS-A
5495307H7G-FO-1	5495307H8G-OOHBA	5495307H9G-SSHB1
5495307H7G-FO-A	5495307H8G-SO-1	5495307H9G-SSHBA
5495307H7G-HH-1	5495307H8G-SO-A	5495307HOG-OOHB1
5495307H7G-HH-A	5495307H8G-SOHB1	5495307HOG-SOHB1
5495307H7G-OO-1	5495307H8G-SOHBA	5497307H1-OO
5495307H7G-OO-A	5495307H8G-SS-1	5497307H1-OOIT
5495307H7G-OOHB1	5495307H8G-SS-A	5497307H1-SO
5495307H7G-OOHBA	5495307H8G-SSHB1	5497307H1-SOIT
5495307H7G-SO-1	5495307H8G-SSHBA	5497307H1-SS
5495307H7G-SO-A	5495307H9-FF-A	5497307H1-SSIT
5495307H7G-SOHB1	5495307H9-OO-A	5497307H1G-OO
5495307H7G-SOHBA	5495307H9-OOHBA	5497307H1G-OOHB

5497307H1G-OOIT	5497307H2G-SS	5498307H1G-SO-1
5497307H1G-SO	5497307H2G-SSHB	5498307H1G-SOHB
5497307H1G-SO-1	5497307H2G-SSHB1	5498307H1G-SOHB1
5497307H1G-SOHB	5497307H2G-SSIT	5498307H1G-SS
5497307H1G-SOHB1	5497307H3-OO	5498307H1G-SSHB
5497307H1G-SOIT	5497307H3-OOIT	5498307H1G-SSHB1
5497307H1G-SOIT1	5497307H3-SO	5498307H2G-OO
5497307H1G-SS	5497307H3-SOIT	5498307H2G-OOHB
5497307H1G-SSHB	5497307H3-SS	5498307H2G-SO
5497307H1G-SSHB1	5497307H3-SSIT	5498307H2G-SO-1
5497307H1G-SSIT	5497307H3G-OO	5498307H2G-SOHB
5497307H2-OO	5497307H3G-OOHB	5498307H2G-SOHB1
5497307H2-OOIT	5497307H3G-OOIT	5498307H2G-SS
5497307H2-SO	5497307H3G-SO	5498307H2G-SS-1
5497307H2-SOIT	5497307H3G-SO-1	5498307H2G-SSHB
5497307H2-SS	5497307H3G-SOHB	5498307H2G-SSHB1
5497307H2-SSIT	5497307H3G-SOHB1	5498307H3G-OO
5497307H2G-OO	5497307H3G-SOIT	5498307H3G-OO-1
5497307H2G-OOHB	5497307H3G-SOIT1	5498307H3G-OOHB
5497307H2G-OOIT	5497307H3G-SS	5498307H3G-SO
5497307H2G-SO	5497307H3G-SSHB	5498307H3G-SO-1
5497307H2G-SO-1	5497307H3G-SSHB1	5498307H3G-SOHB
5497307H2G-SOHB	5497307H3G-SSIT	5498307H3G-SOHB1
5497307H2G-SOHB1	5498307H1G-OO	5498307H3G-SS
5497307H2G-SOIT	5498307H1G-OOHB	5498307H3G-SS-1
5497307H2G-SOIT1	5498307H1G-SO	5498307H3G-SSHB

5498307H3G-SSHB1	5601505G9--B	5605304E0-SO-V
5498307H8-SS-A	5602504E1	5605304E0-SS-V
54B2507G2	5602504E2	5605304E7-HH-A
54B2507G3	5602504E3	5605304E7-OO-1
54B2507G3--G	5602504E3--G	5605304E7-OO-A
54B2807G2	5602505G1	5605304E7-ODPA
54B2807G3	5602505G2	5605304E7-ODPB
54B2807G3--G	5602505G2--G	5605304E7-SO-1
5521607P25	5602505G3	5605304E7-SO-A
5521607P25--B	5602505G3--G	5605304E7-SS-A
5521607P32--B	5602506F1	5605304E8-OO-1
5521607P32-DGB	5602506F3	5605304E8-OO-A
5521607P40	5602604E1	5605304E8-SO-1
5521607P40--B	5602605G3	5605304E8-SO-A
5591207P9--GAK	5602704E1	5605304E8-SS-A
5591507P9--GAB	5602704E3	5605304E9-OO-1
5592507P1--GA	5602705G1	5605304E9-OO-A
5592507P2--GA	5602705G2	5605304E9-SO-1
5592507P3--GA	5602706G3	5605304E9-SO-A
5592707P2--GA	5602804E1	5605304E9-SS-A
5592707P3--GA	5602805G1	5605305G0-OO-1
5592807P1--GA	5602805G2	5605305G0-OO-V
5592807P2--GA	5602805G3	5605305G0-SO-1
5592807P3--GA	5605304E0-OO-1	5605305G0-SO-V
5601504E7--B	5605304E0-OO-V	5605305G0-SS-V
5601504E9--B	5605304E0-SO-1	5605305G7-FF-A

5605305G7-OO-1	5607304E2-SS	5608305G1-OO
5605305G7-OO-A	5607304E3-BB	5608305G1-SO
5605305G7-SO-1	5607304E3-OO	5608305G1-SS
5605305G7-SO-A	5607304E3-SB	5608305G2-OO
5605305G7-SS-A	5607304E3-SO	5608305G2-SO
5605305G8-OO-1	5607304E3-SS	5608305G2-SS
5605305G8-OO-A	5607305G1-OO	5608305G3-OO
5605305G8-SO-1	5607305G1-SO	5608305G3-SO
5605305G8-SO-A	5607305G1-SS	5608305G3-SS
5605305G8-SS-A	5607305G2-FF	5621504E9--B
5605305G8-SS-K	5607305G2-OO	5621505G9--B
5605305G9-OO-1	5607305G2-SO	5621506A9--B
5605305G9-OO-A	5607305G2-SS	5621507A9--B
5605305G9-SO-1	5607305G2-SSDP	5622104E1
5605305G9-SO-A	5607305G3-OO	5622106A1
5605305G9-SS-A	5607305G3-SO	5622106A2
5607304E1-BB	5607305G3-SS	5622107A1
5607304E1-OO	5608304E1-OO	5622107A2
5607304E1-SB	5608304E1-SO	5622107A3
5607304E1-SO	5608304E1-SS	5622503E2
5607304E1-SS	5608304E2-OO	5622503E3
5607304E2-BB	5608304E2-SO	5622503E4
5607304E2-FF	5608304E2-SS	5622504E1
5607304E2-OO	5608304E3-OO	5622504E1--H
5607304E2-SB	5608304E3-SO	5622504E2
5607304E2-SO	5608304E3-SS	5622504E3

5622504E3--C	5622704E1	5625304E7-OO-1
5622504E3--SC	5622704E3	5625304E7-OO-A
5622504E4	5622704E3--C	5625304E7-SO-1
5622504E4--SC	5622705G1	5625304E7-SO-A
5622505G1	5622705G2	5625304E7-SS-A
5622505G2	5622705G3	5625304E7-WW-A
5622505G3	5622706A1	5625304E8-FF-A
5622505G3--SC	5622706F1	5625304E8-OO-1
5622506A1	5622706F3	5625304E8-OO-A
5622506A2	5622707A1	5625304E8-SO-1
5622506A3	5622707A2	5625304E8-SO-A
5622506A3--G	5622707A3	5625304E9-OO-1
5622506F1	5622804E1	5625304E9-OO-A
5622506F2	5622804E2	5625304E9-SO-1
5622506F3	5622804E3	5625304E9-SO-A
5622507A1	5622805G1	5625304E9-SS
5622507A2	5622805G2	5625304E9-SS-A
5622507A3	5622805G3	5625304E9-SS-V
5622603E3	5622806A1	5625305G0-OO-1
5622604E1	5625304E0-OO-1	5625305G0-OO-V
5622604E3	5625304E0-OO-V	5625305G0-SO-1
5622604E3--C	5625304E0-SO-1	5625305G0-SO-V
5622605G3	5625304E0-SO-V	5625305G0-SS-V
5622606F3	5625304E0-SS-V	5625305G7-OO-1
5622606F3--C	5625304E0-SSDPV	5625305G7-SO-1
5622703E3	5625304E4-OO-B	5625305G7-SO-A

5625305G8-OO-1	5625306A9-OO-1	5627304E1-BB
5625305G8-OO-A	5625306A9-OO-A	5627304E1-FF
5625305G8-SO-1	5625306A9-SO-1	5627304E1-OO
5625305G8-SO-A	5625306A9-SS-A	5627304E1-SB
5625305G8-SS-A	5625306F0-OO-V	5627304E1-SO
5625305G9-OO-1	5625306F7-OO-A	5627304E1-SS
5625305G9-OO-A	5625306F8-OO-A	5627304E1-SSDP
5625305G9-SO-1	5625306F9-OO-A	5627304E2-BB
5625305G9-SO-A	5625307A0-FF-V	5627304E2-OO
5625305G9-SS-A	5625307A0-OO-1	5627304E2-SB
5625305G9-WW-A	5625307A0-OO-V	5627304E2-SO
5625306A0-OO-1	5625307A0-SO-1	5627304E2-SS
5625306A0-OO-V	5625307A0-SO-V	5627304E3-BB
5625306A0-SO-1	5625307A0-SS-V	5627304E3-FF
5625306A0-SO-V	5625307A7-OO-1	5627304E3-OO
5625306A0-SS-V	5625307A7-OO-A	5627304E3-SB
5625306A7-FF-A	5625307A7-SO-1	5627304E3-SO
5625306A7-HH-A	5625307A7-SS-A	5627304E3-SS
5625306A7-OO-1	5625307A8-OO-1	5627305G1-BB
5625306A7-OO-A	5625307A8-OO-A	5627305G1-FF
5625306A7-SO-1	5625307A8-SO-1	5627305G1-OO
5625306A7-SS-A	5625307A8-SS-A	5627305G1-SB
5625306A8-OO-1	5625307A9-OO-1	5627305G1-SO
5625306A8-OO-A	5625307A9-OO-A	5627305G1-SS
5625306A8-SO-1	5625307A9-SO-1	5627305G2-BB
5625306A8-SS-A	5625307A9-SS-A	5627305G2-FF

5627305G2-OO	5628305G1-OO	5642503A3--AV
5627305G2-SB	5628305G1-SO	5642503A4--AV
5627305G2-SO	5628305G1-SS	5642703A3--AV
5627305G2-SS	5628305G2-OO	5642803A1--AV
5627305G3-BB	5628305G2-SO	5642803A2--AV
5627305G3-FF	5628305G2-SS	5642803A3
5627305G3-OO	5628305G3-OO	5642803A3--AV
5627305G3-SB	5628305G3-SO	5664002J3--KG
5627305G3-SO	5628305G3-SS	5834005J3--AV
5627305G3-SS	5632504A1--AV	5834005J5--AV
5627306A1-BB	5632504A2--AV	5864003J4--WF
5627306A1-SB	5632504A3--AV	BIOSTAT-KIT1
5627306A2-BB	5632704A3--AV	BIOSTAT-KIT2
5627306A2-SB	5632804A1--AV	BIOSTAT-KIT3
5627306A2-SS	5632804A2--AV	CTR-01
5627306A3-BB	5632804A3--AV	CTR-04
5627306A3-SB	5634002J1	FTR-01
5628304E1-OO	5634002J2	IN-61704-117-K01
5628304E1-SO	5634002J3	IN-61704-119-K01
5628304E1-SS	5634002J4	IN-61704-502-K01
5628304E2-OO	5634003J1	IN-61704-515-O03
5628304E2-SO	5634003J2	IN-61704-517-O03
5628304E2-SS	5634003J3	IN-61704-518-K01
5628304E3-OO	5634003J4	IN-61704-525-O03
5628304E3-SO	5642503A1--AV	IT-CA002PS
5628304E3-SS	5642503A2--AV	IT-CA005GF

J1-5152507T3-E

UNIVESSEL-KIT1

J1-5181307T4-OO

UNIVESSEL-KIT2

J1-5181507T9

J1-5182607T2-LGT

J1-5182658T2-LGT

J1-5185307T8-SS

J1-5185307T9

J1-5185307T9-OO

J1-5231506D9

J1-5231807H9

J1-5235307H7-SS

J1-5235307H9-SS

J1-5441306G5-OO

J1-5441307H4-OO

J1-5445307H7-SS

J1-5445307H8-OO

J1-5601504E9

J1-B65

T5601304E3-SO

T5621305G3-SO

U1-5441307H1OOAC

U1-5441307H4GOOT

U1-5441307H5GOOT

U1-5441358K1SSAC

U1-5441358K2SSAC

U1-5441358K3SSAC
