

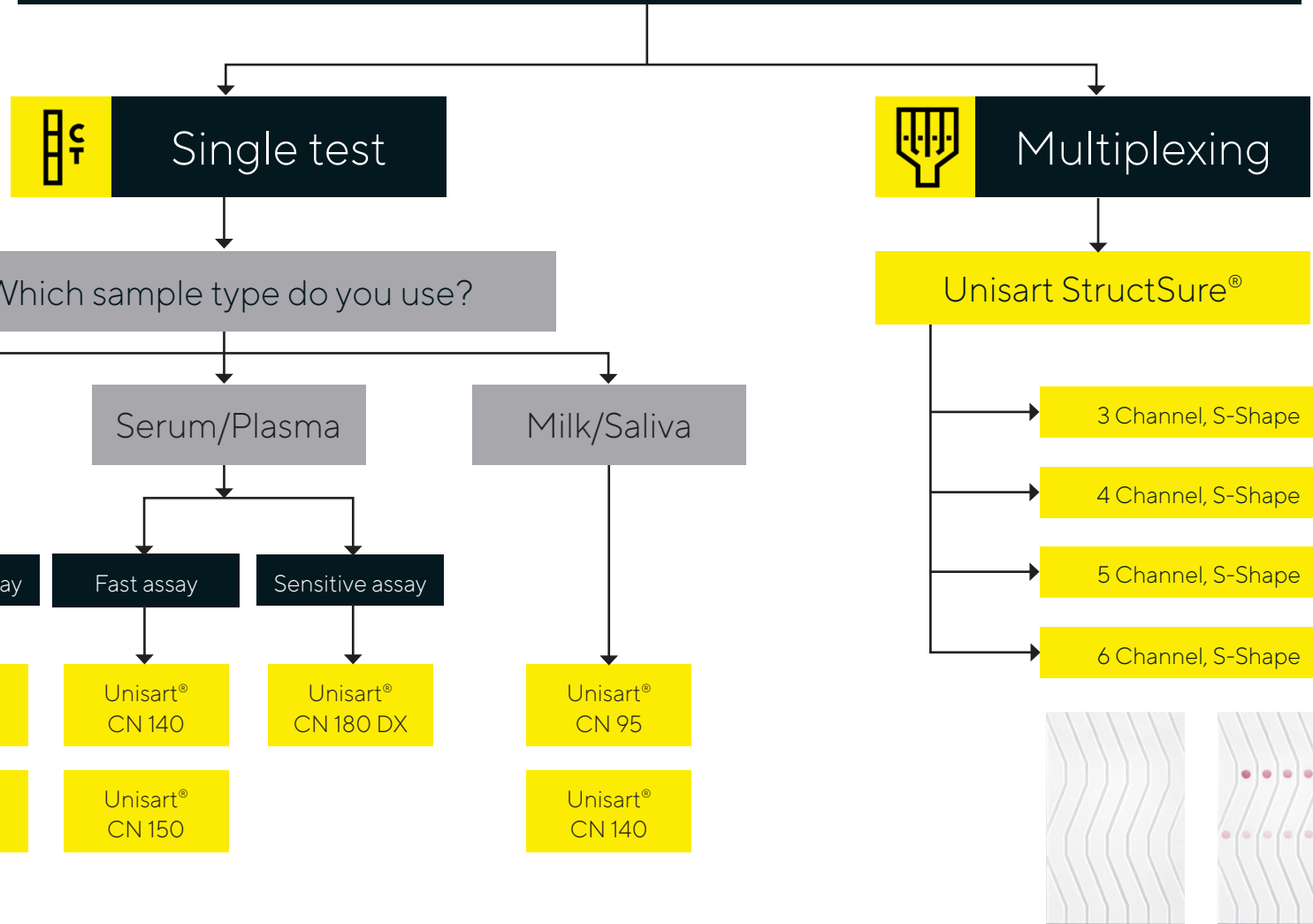


Unisart® CN Membranes  
Selection Guide  
for Lateral Flow Assays

Simplifying Progress

**SARTORIUS**

# Which assay type do you use?



Sartorius offers Unisart® CN Membrane samples for initial evaluation. If you are interested, please contact: [unisart@sartorius.com](mailto:unisart@sartorius.com)

For additional information about Unisart StructSure®, please contact our experts: [unisart@sartorius.com](mailto:unisart@sartorius.com)

All components of a lateral flow test need to be matched and carefully adjusted for every assay, including the heart of every lateral flow test, the membrane. Therefore, it is important to try and select different membrane types according to your individual assay.

Here we provide a suggestion for the membrane with the best fit based on the sample and assay type.

In general, different criteria need to be taken under consideration such as

- Capillary speed
- Sensitivity
- Sample viscosity
- Antibody-antigen system
- Binding kinetics
- Label type
- Reader system required



Please visit our website for more information or ask our experts for support ([unisart@sartorius.com](mailto:unisart@sartorius.com)).



Ready to buy?  
Order Unisart® CN Membranes online via the Sartorius eShop.

## Specifications and Characteristics

Unisart®	CN 95 backed	CN 110 backed	CN 140 unbacked	CN 140 backed	CN 150 backed	CN 180 DX backed
Membrane material	Nitrocellulose	Nitrocellulose	Nitrocellulose	Nitrocellulose	Nitrocellulose	Nitrocellulose
Backing	100 µm polyester, clear	100 µm polyester, clear	none	100 µm polyester, clear	100 µm polyester, clear; or 100 µm polyester, white	100 µm polyester, clear
Thickness (µm)	240 - 270	185 - 215	120 - 160	225 - 255	240 - 280	225 - 255
Capillary speed down web, purified water (s/40mm)	65 - 115	90 - 130	90 - 150	95 - 155	90 - 180	135 - 175

## Ordering Information

Unisart® Membrane	Polyester backing	Width (mm)	Length (m)	Max. splices	Units   Box	Catalog number
<b>CN 95 backed</b>						
CN 95	100 µm clear	20	100	3	1 roll	1UN95ER100020NT
CN 95	100 µm clear	22	100	3	5 rolls	1UN95ER100022NTB
CN 95	100 µm clear	25	100	3	5 rolls	1UN95ER100025NTB
CN 95	100 µm clear	30	100	3	5 rolls	1UN95ER100030-B
CN 95	100 µm clear	35	100	3	5 rolls	1UN95ER100035NTB
CN 95	100 µm clear	40	100	0	1 roll	1UN95ER100040WS
<b>CN 110 backed</b>						
CN 110	100 µm clear	20	100	3	5 rolls	1UN11ER100020NTB
CN 110	100 µm clear	25	100	3	1 roll	1UN11ER100025NT
CN 110	100 µm clear	25	100	3	5 rolls	1UN11ER100025NTB
<b>CN 140 unbacked</b>						
CN 140	none	20	100	3	5 rolls	1UN14AR100020NTB
CN 140	none	20	100	0	5 rolls	1UN14AR100020WSB
CN 140	none	25	100	3	5 rolls	1UN14AR100025NTB
CN 140	none	25	100	0	5 rolls	1UN14AR100025WSB
CN 140	none	35	100	3	5 rolls	1UN14AR100035NTB
CN 140	none	270	100	3	1 roll	1UN14AR100270NT
<b>CN 140 backed</b>						
CN 140	100 µm clear	18	100	3	5 rolls	1UN14ER100018NTB
CN 140	100 µm clear	20	100	3	1 roll	1UN14ER100020NT
CN 140	100 µm clear	20	100	3	5 rolls	1UN14ER100020NTB
CN 140	100 µm clear	25	100	3	1 roll	1UN14ER100025NT
CN 140	100 µm clear	25	100	3	5 rolls	1UN14ER100025NTB
CN 140	100 µm clear	35	100	3	5 rolls	1UN14ER100035NTB
CN 140	100 µm clear	40	100	0	5 rolls	1UN14ER100040WSB

Customized dimensions of Unisart® CN Membranes are also available upon request.

For further information and customized dimensions please contact us: [unisart@sartorius.com](mailto:unisart@sartorius.com)

Some products can be ordered online via the [Sartorius eShop](#).

Unisart® Membrane	Polyester backing	Width (mm)	Length (m)	Max. splices	Units   Box	Catalog number
<b>CN 150 backed</b>						
CN 150	100 µm clear	20	100	3	5 rolls	1UN15LR100020NTB
CN 150	100 µm clear	25	100	3	1 roll	1UN15LR100025NT
CN 150	100 µm clear	25	100	3	5 rolls	1UN15LR100025NTB
CN 150	100 µm white	25	100	3	1 roll	1UN15WR100025NT
CN 150	100 µm white	25	100	3	5 rolls	1UN15WR100025NTB
<b>CN 180 DX backed</b>						
CN 180 DX	100 µm clear	25	100	3	1 roll	1DX18ER100020NT
CN 180 DX	100 µm clear	20	100	3	5 rolls	1DX18ER100020NTB
CN 180 DX	100 µm clear	25	100	3	1 roll	1DX18ER100025NT
CN 180 DX	100 µm clear	25	100	3	5 rolls	1DX18ER100025NTB
<b>StructSure®</b>						
3 Channel, S-Shape	100 µm clear	25	100	0	1 roll	3UN14ER103S01WS
3 Channel, S-Shape	100 µm clear	25	100	0	5 rolls	3UN14ER103S01WSB
4 Channel, S-Shape	100 µm clear	25	100	0	1 roll	3UN14ER084S01WS
4 Channel, S-Shape	100 µm clear	25	100	0	5 rolls	3UN14ER084S01WSB
5 Channel, S-Shape	100 µm clear	25	100	0	1 roll	3UN14ER075S01WS
5 Channel, S-Shape	100 µm clear	25	100	0	5 rolls	3UN14ER075S01WSB
6 Channel, S-Shape	100 µm clear	25	100	0	1 roll	3UN14ER066S01WS
6 Channel, S-Shape	100 µm clear	25	100	0	5 rolls	3UN14ER066S01WSB

Customized dimensions of Unisart® CN Membranes are also available upon request.  
For further information and customized dimensions please contact us: [unisart@sartorius.com](mailto:unisart@sartorius.com)  
Some products can be ordered online via the [Sartorius eShop](#).



## Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Or +1 650 322 1360



For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)