according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

SECTION 1: Identification of the substance/mixture and of the company/undertaking

1.1 Product identifier

Trade name : 4Cell MDCK CD Medium 10L Powder

Product code : CQV3FA2012

Substance name : 4Cell® MDCK CD Powder Medium

1.2 Relevant identified uses of the substance or mixture and uses advised against

1.3 Details of the supplier of the safety data sheet

Company : Sartorius Stedim Biotech GmbH

August-Spindler-Strasse 11

37079 Göttingen

E-mail address of person

responsible for the SDS

: Department CA-FM Environment@sartorius.com

1.4 Emergency telephone number

UK/Europe/Global: +44.1865.407.333 (Sartorius 29003)

United States: +1.866.928.0789 (Toll free), (Sartorius 29003)

Canada: +1.800.579.7421 (Toll free), (Sartorius 29003)

United States & Canada (geographic): +1.202.464.2554 (Sartorius 29003)

Asia-Pacific/East & South East Asia (without China): +65.3158.1074 (Sartorius 29003)

Japan: +81.3.4578.9341 (Sartorius 29003)

Indonesia: 007.803.011.0293 (Toll free Indonesia), (Sartorius 29003)

Malaysia: +60.3.6207.4347 (Sartorius 29003)

Thailand: 001.800.120.666.751 (Toll free Thailand), (Sartorius 29003)

India: 000.800.100.7479 (Toll free India) or +65.3158.1198 (Sartorius 29003)

South Korea: +82.2.3479.8401 (Sartorius 29003) Australia: +61.2.8014.4558 (Sartorius 29003)

New Zealand: +64.9.929.1483 (Sartorius 29003)

Middle East/Africa: +44.1235.239671 (Sartorius 29003)

SECTION 2: Hazards identification

2.1 Classification of the substance or mixture

Classification (REGULATION (EC) No 1272/2008)

Eye irritation, Category 2 H319: Causes serious eye irritation.

Skin irritation, Category 2 H315: Causes skin irritation.

Specific target organ toxicity - single ex- H335: May cause respiratory irritation.

posure, Category 3

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

2.2 Label elements

Labelling REGULATION (EC) No 1272/2008

Hazard pictograms

Signal word : Warning

Hazard statements : H319 Causes serious eye irritation.

H315 Causes skin irritation.

H335 May cause respiratory irritation.

Precautionary statements : Prevention:

P261 Avoid breathing dust/ fume/ gas/ mist/ va-

pours/ spray.

P271 Use only outdoors or in a well-ventilated

area.

P280 Wear protective gloves/ protective clothing/

eye protection/ face protection.

Response:

P305 + P351 + P338 IF IN EYES: Rinse cautiously with wa-

ter for several minutes. Remove contact lenses, if present and easy to do. Continue

rinsing.

P337 + P313 If eye irritation persists: Get medical advice/

attention.

P302 + P352 IF ON SKIN: Wash with plenty of soap and

water.

P332 + P313 If skin irritation occurs: Get medical advice/

attention.

P304 + P340 IF INHALED: Remove victim to fresh air

and keep at rest in a position comfortable

for breathing.

P312 Call a POISON CENTER/doctor if you feel

unwell.

Storage:

P403 + P233 Store in a well-ventilated place. Keep con-

tainer tightly closed.

2.3 Other hazards

None known.

SECTION 3: Composition/information on ingredients

3.2 Mixtures

Chemical nature : Reagent

Mixture

powdered form

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

Hazardous components

CAS-No. EC-No. Registration number	Chemical name	Classification (67/548/EEC)	Classification (REGULATION (EC) No 1272/2008)	Concentration (%)
10102-18-8	Sodium sele- nite 233-267-9	T; T; R23/25 N; R51 R31	Acute Tox. 2; H300 Acute Tox. 2; H330 Skin Irrit. 2; H315 6.4 2; H319 Skin Sens. 1; H317 Aquatic Chronic 2; H411	<= 0,1
110-15-6	Succinic acid 203-740-4		Eye Dam. 1; H318	<= 0,1
7446-20-0	Zinc sulfate heptahydrate 616-097-3		Acute Tox. 4; H302 Eye Dam. 1; H318 Aquatic Acute 1; H400 Aquatic Chronic 1; H410	<= 0,1
30925-07-6	L-Cystine dihy- drochloride 250-391-9		Skin Corr. 1B; H314 Eye Dam. 1; H318	<= 1

For explanation of abbreviations see section 16.

SECTION 4: First aid measures

4.1 Description of first aid measures

If inhaled : Move to fresh air.

If symptoms persist, call a physician.

In case of skin contact : Wash off immediately with soap and plenty of water.

In case of eye contact : Rinse immediately with plenty of water, also under the eyelids.

If eye irritation persists, consult a specialist.

If swallowed : Rinse mouth with water.

Never give anything by mouth to an unconscious person.

Consult a physician if necessary.

4.2 Most important symptoms and effects, both acute and delayed

None known.

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019 18.05.2020 12

4.3 Indication of any immediate medical attention and special treatment needed

Treatment : No information available.

SECTION 5: Firefighting measures

5.1 Extinguishing media

Suitable extinguishing media Use water spray, alcohol-resistant foam, dry chemical or car-

bon dioxide.

5.2 Special hazards arising from the substance or mixture

Specific hazards during fire-

fighting

: No information available.

5.3 Advice for firefighters

for firefighters

Special protective equipment : In the event of fire, wear self-contained breathing apparatus.

Further information : The product itself does not burn.

Use extinguishing measures that are appropriate to local cir-

cumstances and the surrounding environment.

SECTION 6: Accidental release measures

6.1 Personal precautions, protective equipment and emergency procedures

Personal precautions : Avoid dust formation.

6.2 Environmental precautions

Environmental precautions : Do not let product enter drains.

6.3 Methods and material for containment and cleaning up

: Pick up and transfer to properly labelled containers. Methods for cleaning up

6.4 Reference to other sections

SECTION 7: Handling and storage

7.1 Precautions for safe handling

Hygiene measures : Avoid contact with skin, eyes and clothing. Do not breathe

dust. Keep away from food, drink and animal feedingstuffs.

When using, do not eat, drink or smoke.

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

7.2 Conditions for safe storage, including any incompatibilities

Requirements for storage areas and containers

: Keep container tightly closed. Keep in the refrigerator (2-8°C).

7.3 Specific end use(s)

SECTION 8: Exposure controls/personal protection

8.1 Control parameters

8.2 Exposure controls

Personal protective equipment

Eye protection

: Goggles

Hand protection

Material : Nitrile rubber

butyl-rubber

Skin and body protection : Lab coat

SECTION 9: Physical and chemical properties

9.1 Information on basic physical and chemical properties

Appearance : powder

Colour : Faint yellow to faint red

Odour : specific smell

pH : ≥ 3,28

to ≤ 3,92

9.2 Other information

No data available

SECTION 10: Stability and reactivity

10.1 Reactivity

10.2 Chemical stability

10.3 Possibility of hazardous reactions

Hazardous reactions : No dangerous reaction known under conditions of normal use.

No decomposition if stored and applied as directed.

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

10.4 Conditions to avoid

10.5 Incompatible materials

Materials to avoid : Oxidizing agents

10.6 Hazardous decomposition products

No decomposition if stored and applied as directed.

SECTION 11: Toxicological information

11.1 Information on toxicological effects

Further information

Product:

Remarks: Health injuries are not known or expected under normal use.

SECTION 12: Ecological information

12.1 Toxicity

No data available

12.2 Persistence and degradability

No data available

12.3 Bioaccumulative potential

No data available

12.4 Mobility in soil

No data available

12.5 Results of PBT and vPvB assessment

Not relevant

12.6 Other adverse effects

Product:

Additional ecological infor-

: Remarks: This product has no known ecotoxicological effects.

mation

SECTION 13: Disposal considerations

13.1 Waste treatment methods

Product : In accordance with local and national regulations.

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

SECTION 14: Transport information

- 14.1 UN number
- 14.2 UN proper shipping name
- 14.3 Transport hazard class(es)
- 14.4 Packing group
- 14.5 Environmental hazards
- 14.6 Special precautions for user

Not applicable

14.7 Transport in bulk according to Annex II of MARPOL 73/78 and the IBC Code

Not applicable for product as supplied.

SECTION 15: Regulatory information

15.1 Safety, health and environmental regulations/legislation specific for the substance or mixture

Water contaminating class

(Germany)

: WGK 1 slightly hazardous to water

15.2 Chemical safety assessment SECTION 16: Other information

Full text of R-Phrases

R23/25 : Toxic by inhalation and if swallowed.
R31 : Contact with acids liberates toxic gas.

R51 : Toxic to aquatic organisms.

Full text of H-Statements

H300 : Fatal if swallowed. H302 : Harmful if swallowed.

H314 : Causes severe skin burns and eye damage.

H315 : Causes skin irritation.

H317 : May cause an allergic skin reaction.
H318 : Causes serious eye damage.
H319 : Causes serious eye irritation.

H330 : Fatal if inhaled.

H400 : Very toxic to aquatic life.

H410 : Very toxic to aquatic life with long lasting effects.H411 : Toxic to aquatic life with long lasting effects.

Full text of other abbreviations

according to Regulation (EC) No. 1907/2006

4Cell MDCK CD Medium 10L Powder

Version Revision Date: SDS Number: Date of last issue: 18.05.2020 1.2 18.05.2020 SAG-R0002081 Date of first issue: 16.12.2019

6.4 : Eye irritation Acute Tox. : Acute toxicity

Aquatic Acute : Short-term (acute) aquatic hazard Aquatic Chronic : Long-term (chronic) aquatic hazard

Eye Dam. : Serious eye damage
Skin Corr. : Skin corrosion
Skin Irrit. : Skin irritation
Skin Sens. : Skin sensitisation

The information provided in this Safety Data Sheet is correct to the best of our knowledge, information and belief at the date of its publication. The information given is designed only as a guidance for safe handling, use, processing, storage, transportation, disposal and release and is not to be considered a warranty or quality specification. The information relates only to the specific material designated and may not be valid for such material used in combination with any other materials or in any process, unless specified in the text.

DE / EN