

SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	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 Liquid Medium
Substance name : 4Cell® MDCK CD Liquid 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)

Skin irritation, Category 2	H316: Causes mild skin irritation.
Eye irritation, Category 2	H320: Causes eye irritation.
Specific target organ toxicity - single exposure, Category 3	H335: May cause respiratory irritation.

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	Date of first issue: 16.12.2019

2.2 Label elements

Labelling REGULATION (EC) No 1272/2008

Hazard pictograms :



Signal word : Warning

Hazard statements : H316 Causes mild skin irritation.
H320 Causes eye irritation.
H335 May cause respiratory irritation.

Precautionary statements : P261 Avoid breathing dust/ fume/ gas/ mist/ va-
pours/ spray.
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 + P341 IF INHALED: If breathing is difficult, remove
victim to fresh air and keep at rest in a posi-
tion comfortable for breathing.
P312 Call a POISON CENTER/doctor/.? if you
feel unwell.

2.3 Other hazards

EUH208: Contains sodium selenite. May cause allergic reactions.

SECTION 3: Composition/information on ingredients

3.2 Mixtures

Chemical nature : Aqueous solution
Reagent

Hazardous components

CAS-No. EC-No. Registration number	Chemical name	Classification (67/548/EEC)	Classification (REGULATION (EC) No	Concentration (%)
--	---------------	--------------------------------	--	----------------------

4Cell® MDCK CD Liquid Medium

Version 1.3 Revision Date: 18.05.2020 SDS Number: SAG-R0002082 Date of last issue: 09.03.2020
Date of first issue: 16.12.2019

			1272/2008)	
1310-73-2	Sodium hydroxide 215-185-5	C-Xi; C-Xi; R35-R41	Met. Corr. 1; H290 Skin Corr. 1A; H314 Eye Dam. 1; H318	<= 0,01
10102-18-8	Sodium selenite 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	<= 0,01
110-15-6	Succinic acid 203-740-4		Eye Dam. 1; H318	<= 0,01
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	<= 0,01
30925-07-6	L-Cystine dihydrochloride 250-391-9		Skin Corr. 1B; H314 Eye Dam. 1; H318	<= 0,1

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.

4.3 Indication of any immediate medical attention and special treatment needed

Treatment : No information available.

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	Date of first issue: 16.12.2019

SECTION 5: Firefighting measures

5.1 Extinguishing media

Suitable extinguishing media : Use water spray, alcohol-resistant foam, dry chemical or carbon dioxide.

5.2 Special hazards arising from the substance or mixture

Specific hazards during fire-fighting : No information available.

5.3 Advice for firefighters

Special protective equipment for firefighters : 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 circumstances and the surrounding environment.

SECTION 6: Accidental release measures

6.1 Personal precautions, protective equipment and emergency procedures

Personal precautions : Avoid formation of aerosol.

6.2 Environmental precautions

Environmental precautions : Do not let product enter drains.

6.3 Methods and material for containment and cleaning up

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

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. Keep away from food, drink and animal feedingstuffs. When using do not eat or drink.

7.2 Conditions for safe storage, including any incompatibilities

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

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	Date of first issue: 16.12.2019

areas and containers : Protect from light.

Storage class (TRGS 510) : 10-13, German Storage Class 10 to 13

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	: butyl-rubber Nitrile rubber
Skin and body protection	: Lab coat

SECTION 9: Physical and chemical properties

9.1 Information on basic physical and chemical properties

Appearance	: liquid, Aqueous solution
Colour	: clear yellow to slightly red
Odour	: specific smell
pH	: $\geq 7,0$ to $\leq 7,3$
Density	: ca. 1 g/cm ³ similar to water

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.

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	Date of first issue: 16.12.2019

No decomposition if stored and applied as directed.

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 information : Remarks: In large quantities, this product may affect the functioning of waste-water treatment plants.

SECTION 13: Disposal considerations

13.1 Waste treatment methods

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	Date of first issue: 16.12.2019

Product : In accordance with local and national regulations.

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

Remarks : Not dangerous goods in the meaning of ADR/RID, ADN, IMDG-Code, ICAO/IATA-DGR

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.
R35	: Causes severe burns.
R41	: Risk of serious damage to eyes.
R51	: Toxic to aquatic organisms.

Full text of H-Statements

H290	: May be corrosive to metals.
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.

SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

4Cell® MDCK CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 09.03.2020
1.3	18.05.2020	SAG-R0002082	Date of first issue: 16.12.2019

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

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
Met. Corr.	: Corrosive to metals
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