

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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 100L Powder

Product code : CQV3FA2014

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 exposure, Category 3	H335: May cause respiratory irritation.

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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/ vapours/ 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 water 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 container tightly closed.

### 2.3 Other hazards

None known.

---

## SECTION 3: Composition/information on ingredients

### 3.2 Mixtures

Chemical nature : Reagent  
Mixture  
powdered form

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

## 4Cell MDCK CD Medium 100L Powder

Version 1.2      Revision Date: 18.05.2020      SDS Number: SAG-R0002081      Date of last issue: 18.05.2020  
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 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	<= 0,1
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 dihydrochloride 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.

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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

---

### 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 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 dust formation.

### 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. Do not breathe dust. Keep away from food, drink and animal feedingstuffs. When using, do not eat, drink or smoke.

---

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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	: $\geq 3,28$  to $\leq 3,92$

### 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.

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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 information : Remarks: This product has no known ecotoxicological effects.

---

## SECTION 13: Disposal considerations

### 13.1 Waste treatment methods

Product : In accordance with local and national regulations.

---

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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 : WGK 1 slightly hazardous to water  
(Germany)

#### 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

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

---

## 4Cell MDCK CD Medium 100L 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