

# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

---

## 4Cell® BHK-21 CD Powder Medium

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

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

#### 1.1 Product identifier

Trade name : 4Cell® BHK-21 CD Powder Medium  
Substance name : 4Cell® BHK-21 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)

Skin irritation, Category 2 H315: Causes skin irritation.

Eye irritation, Category 2 H320: Causes eye irritation.

Specific target organ toxicity - single exposure, Category 2 H335: May cause respiratory irritation.

---


## 4Cell® BHK-21 CD Powder Medium

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

### 2.2 Label elements

#### Labelling REGULATION (EC) No 1272/2008

Hazard pictograms	:		
Signal word	:	Warning	
Hazard statements	:	H315 H320 H335	Causes skin irritation. Causes eye irritation. May cause respiratory irritation.
Precautionary statements	:	<b>Prevention:</b> P261 P280 <b>Response:</b> P305 + P351 + P338 P337 + P313 P302 + P352 P332 + P313 P304 + P341 P312	Avoid breathing dust. Wear protective gloves/ protective clothing/ eye protection/ face protection. IF IN EYES: Rinse cautiously with wa- ter for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. If eye irritation persists: Get medical advice/ attention. IF ON SKIN: Wash with plenty of soap and water. If skin irritation occurs: Get medical advice/ attention. IF INHALED: If breathing is difficult, remove victim to fresh air and keep at rest in a posi- tion comfortable for breathing. Call a POISON CENTER/doctor/.? if you feel unwell.

### 2.3 Other hazards

EUH208: Contains cobalt (II) chloride hexahydrate, sodium pyruvate and sodium selenite. Can cause allergic reactions.

## SECTION 3: Composition/information on ingredients

### 3.2 Mixtures

Chemical nature	:	Reagent powdered form
-----------------	---	--------------------------

#### Hazardous components

**SAFETY DATA SHEET**

according to Regulation (EC) No. 1907/2006

**SARTORIUS****4Cell® BHK-21 CD Powder Medium**Version  
1.3Revision Date:  
15.04.2020SDS Number:  
SAG-R0001802Date of last issue: 21.02.2020  
Date of first issue: 05.03.2019

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
7791-13-1	Cobalt chloride hexahydrate	Xn-N; Xn; R45- R20/21/22-R42- R52/53	Acute Tox. 4; H302 Skin Sens. 1; H317 Resp. Sens. 1; H334 Muta. 2; H341 Carc. 1B; H350i Repr. 1B; H360F Aquatic Acute 1; H400 Aquatic Chronic 1; H410	<= 0,1
7803-55-6	Ammonium metavanadate 232-261-3	T; T; R25 T; T; R48-R23 Repr.Cat.3; R62- R63 Xn; Xn; R20 N; R51/53	Acute Tox. 3; H301 Acute Tox. 4; H332 Eye Irrit. 2; H319 Repr. 2; H361 STOT RE 1; H372 Aquatic Chronic 2; H411	<= 0,1
10025-69-1	Tin chloride dihydrate	Xn; Xn; R21 C; C; R34 N; N; R50/53	Acute Tox. 4; H302 Skin Corr. 1B; H314 Aquatic Acute 1; H400	<= 0,1
13517-24-3	Sodium meta- silicate no- nahydrate 229-912-9		Met. Corr. 1; H290 Skin Corr. 1B; H314 Eye Dam. 1; H318 STOT SE 3; H335	<= 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
113-24-6	Sodium py- ruvate 204-024-4		Eye Irrit. 2; H319 Skin Sens. 1B; H317	<= 0,9
45738-97-4	3,7-		Eye Dam. 1; H318	<= 0,1

---

## **4Cell® BHK-21 CD Powder Medium**

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

	dihydropurin-6-one, Sodium			
--	----------------------------	--	--	--

For explanation of abbreviations see section 16.

### **SECTION 4: First aid measures**

#### **4.1 Description of first aid measures**

- If inhaled : If breathed in, move person into fresh air.  
In the case of inhalation of aerosol/mist consult a physician if necessary.
- In case of skin contact : Wash off immediately with soap and plenty of water.  
If skin irritation persists, call a physician.
- In case of eye contact : In the case of contact with eyes, rinse immediately with plenty of water and seek medical advice.
- If swallowed : Clean mouth with water and drink afterwards plenty of water.  
Never give anything by mouth to an unconscious person.  
Call a physician immediately.

#### **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**

### **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**

#### **5.3 Advice for firefighters**

- Special protective equipment for firefighters : Wear self-contained breathing apparatus for firefighting if necessary.
- Further information : The product itself does not burn.  
Use extinguishing measures that are appropriate to local circumstances and the surrounding environment.

---

## **4Cell® BHK-21 CD Powder Medium**

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

### **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 arrange disposal without creating dust.  
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. Wear suitable gloves and eye/face protection.

#### **7.2 Conditions for safe storage, including any incompatibilities**

Requirements for storage areas and containers : Keep container tightly closed in a dry and well-ventilated place. Keep in the refrigerator (2-8°C).

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 : Tightly fitting safety goggles

Hand protection

Remarks : butyl-rubber Nitrile rubber

Skin and body protection : Lab coat

---

## **4Cell® BHK-21 CD Powder Medium**

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

### **SECTION 9: Physical and chemical properties**

#### **9.1 Information on basic physical and chemical properties**

Appearance : solid

Colour : colourless, white, reddish-violet

Odour : odourless

Density : No data available

#### **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.  
Stable under recommended storage conditions.

#### **10.4 Conditions to avoid**

Conditions to avoid : Exposure to moisture

#### **10.5 Incompatible materials**

Materials to avoid : Strong oxidizing agents

#### **10.6 Hazardous decomposition products**

### **SECTION 11: Toxicological information**

#### **11.1 Information on toxicological effects**

##### **Further information**

##### **Product:**

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

---

## **4Cell® BHK-21 CD Powder Medium**

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

### **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**

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.

---

## **4Cell® BHK-21 CD Powder Medium**

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

### **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**

R20	: Harmful by inhalation.
R20/21/22	: Harmful by inhalation, in contact with skin and if swallowed.
R21	: Also harmful in contact with skin.
R23	: Toxic by inhalation.
R23/25	: Toxic by inhalation and if swallowed.
R25	: Toxic if swallowed.
R31	: Contact with acids liberates toxic gas.
R34	: Causes burns.
R42	: May cause sensitisation by inhalation.
R45	: May cause cancer.
R48	: Danger of serious damage to health by prolonged exposure.
R50/53	: Very toxic to aquatic organisms, may cause long-term adverse effects in the aquatic environment.
R51	: Toxic to aquatic organisms.
R51/53	: Toxic to aquatic organisms, may cause long-term adverse effects in the aquatic environment.
R52/53	: Harmful to aquatic organisms, may cause long-term adverse effects in the aquatic environment.
R62	: Possible risk of impaired fertility.
R63	: Possible risk of harm to the unborn child.

#### **Full text of H-Statements**

H290	: May be corrosive to metals.
H300	: Fatal if swallowed.
H301	: Toxic 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.
H332	: Harmful if inhaled.
H334	: May cause allergy or asthma symptoms or breathing difficulties if inhaled.



# SAFETY DATA SHEET

according to Regulation (EC) No. 1907/2006

SARTORIUS

---

## 4Cell® BHK-21 CD Powder Medium

Version	Revision Date:	SDS Number:	Date of last issue: 21.02.2020
1.3	15.04.2020	SAG-R0001802	Date of first issue: 05.03.2019

---

H335	: May cause respiratory irritation.
H341	: Suspected of causing genetic defects if inhaled.
H350i	: May cause cancer by inhalation.
H360F	: May damage fertility.
H361	: Suspected of damaging fertility or the unborn child.
H372	: Causes damage to organs through prolonged or repeated exposure 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
Carc.	: Carcinogenicity
Eye Dam.	: Serious eye damage
Eye Irrit.	: Eye irritation
Met. Corr.	: Corrosive to metals
Muta.	: Germ cell mutagenicity
Repr.	: Reproductive toxicity
Resp. Sens.	: Respiratory sensitisation
Skin Corr.	: Skin corrosion
Skin Irrit.	: Skin irritation
Skin Sens.	: Skin sensitisation
STOT RE	: Specific target organ toxicity - repeated exposure
STOT SE	: Specific target organ toxicity - single exposure

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