SVIFCTFAS

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

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

1.1 Product identifier

Trade name	:	4Cell® BHK-21 CD Liquid Medium
Substance name	:	4Cell® BHK-21 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.
Skin irritation, Category 2	H320: Causes eye irritation.
Specific target organ toxicity - single ex- posure, Category 2	H335: May cause respiratory irritation.

SVILCILVS

according to Regulation (EC) No. 1907/2006

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

2.2 Label elements

Labelling REGULATION (EC)	No 1272/2008	
Hazard pictograms :		
Signal word :	Warning	
Hazard statements :	H316 H320 H335	Causes mild skin irritation. Causes eye irritation. May cause respiratory irritation.
Precautionary statements :	P261 P280	Avoid breathing spray. Wear protective gloves/ protective clothing/ eye protection/ face protection.
	Response: P305 + P351 + P3	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 position comfortable for breathing.
	P312	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

: Mixture Aqueous solution

Hazardous components

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

SVILOTEVS

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date: 14.12.2020	SDS Number:	Date of last issue: 30.04.2020
1.1		SAG-R0002080	Date of first issue: 30.04.2020

			1272/2008)	
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	<= 0,01
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	<= 0,01
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	<= 0,01
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	<= 0,01
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	<= 0,01
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
113-24-6	Sodium py- ruvate 204-024-4		Eye Irrit. 2; H319 Skin Sens. 1B; H317	<= 0,1
45738-97-4	3,7- dihydropurin-6- one, Sodium		Eye Dam. 1; H318	<= 0,01

SVIFCTSVS

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

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 thoroughly 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.

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- : No information available. fighting

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 cir- cumstances and the surrounding environment.	

SVIFCTFAS

according to Regulation (EC) No. 1907/2006

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

SECTION 6: Accidental release measures

6.1 Personal precautions, protective equipment and emergency procedures			
Personal precautions	: Avoid formation of aerosol.		
6.2 Environmental precautions			
•	: Should not be released into the environment.		
Environmental precautions	Do not let product enter drains.		
6.3 Methods and material for con	tainment 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			
SECTION 7. Handling and Storage			

7.1 Precautions for safe handling Advice on safe handling Hygiene measures Avoid formation of aerosol. Avoid contact with skin, eyes and clothing. Keep away from food and drink. When using do not eat or drink.

7.2 Conditions for safe storage, including any incompatibilities Requirements for storage areas and containers Protect from exposure to light. Keep container tightly closed. 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

No data available

8.2 Exposure controls

Personal protective equipment		
Eye protection :	Goggles	
Hand protection		
Remarks :	butyl-rubber Nitrile rubber	
Skin and body protection :	Lab coat	

SVILCIEVS

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

SECTION 9: Physical and chemical properties

9.1 Information on basic physical and chemical properties

Appearance Colour Odour pH	 liquid, Aqueous solution clear, light yellow specific smell 6,9
F	to ≤ 7,1
Density	: ca. 1 g/cm3 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	:	None known.
		No decomposition if stored and applied as directed.

10.4 Conditions to avoid

10.5 Incompatible materials

Materials to avoid

: Oxidizing agents Acids and bases

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.

SVIFCTFAS

according to Regulation (EC) No. 1907/2006

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

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: In large quantities, this product may affect the func-
mation		tioning 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.

SVIECTENS

according to Regulation (EC) No. 1907/2006

4Cell® BHK-21 CD Liquid Medium

Version	Revision Date:	SDS Number:	Date of last issue: 30.04.2020
1.1	14.12.2020	SAG-R0002080	Date of first issue: 30.04.2020

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 R20/21/22 R21 R23 R23/25 R25 R31 R34 R42 R45 R48 R50/53 R51 R51/53 R52/53	 Harmful by inhalation. Harmful by inhalation, in contact with skin and if swallowed. Also harmful in contact with skin. Toxic by inhalation. Toxic by inhalation and if swallowed. Toxic if swallowed. Contact with acids liberates toxic gas. Causes burns. May cause sensitisation by inhalation. May cause cancer. Danger of serious damage to health by prolonged exposure. Very toxic to aquatic organisms, may cause long-term adverse effects in the aquatic environment. Toxic to aquatic organisms, may cause long-term adverse effects in the aquatic environment. Harmful to aquatic organisms, may cause long-term adverse effects in the aquatic environment.
R62 R63	Possible risk of impaired fertility.Possible risk of harm to the unborn child.
Full text of H-Statements	
H290	: May be corrosive to metals.
H300 H301	: Fatal if swallowed. : 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 difficul- ties if inhaled.

SVILOTEVS

4Cell® BHK-21 CD Liquid Medium

Version 1.1	Revision Date: 14.12.2020	SDS Number: SAG-R0002080	Date of last issue: 30.04.2020 Date of first issue: 30.04.2020
H335 H341 H350i H360F H361 H372 H400 H410	-	 Suspected of c May cause car May damage f Suspected of c Causes damage exposure if inh Very toxic to an Very toxic to an 	lamaging fertility or the unborn child. ge to organs through prolonged or repeated aled. quatic life. quatic life with long lasting effects.
H411 Full te	ext of other abbreviati	•	c life with long lasting effects.
6.4 Acute Aquati	Tox. c Acute c Chronic am. it. corr. Sens. corr. rit. eens. RE	 Eye irritation Acute toxicity Short-term (ac Long-term (chr Carcinogenicity Serious eye da Eye irritation Corrosive to m Germ cell muta Reproductive t Respiratory se Skin corrosion Skin sensitisat Specific target 	amage etals agenicity oxicity nsitisation

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