

1. Identification of the substance/preparation and the company/undertaking

Identification of the substance/preparation

Product Code: NBP2-31290 thru NBP2-31294

Product Name: Novus Neural Human Progenitor Cells (hPNCs)

Company/Undertaking Identification

Novus Biologicals, LLC
8100 Southpark Way, A-8
Littleton, CO 80120
USA

2. Composition/Information on ingredients

• Hazardous/Non-hazardous Components:

Less than 10% of Dimethyl sulfoxide (DMSO) is present in frozen cultures.

Dimethyl sulfoxide (DMSO)

CAS NO: 67-68-5

MF: C₂H₆SO

EC NO: 200-664-3

Contact Novus Biologicals, LLC for further information. All other ingredients in the proportions used are considered non hazardous by regulating bodies worldwide.

3. Hazards Identification

- **Classification/ Indication of danger:** Not hazardous

4. First Aid Measures

- **Skin Contact:** Wash off immediately with plenty of water.
- **Eye Contact:** Rinse thoroughly with plenty of water, also under eyelids.
- **Ingestion:** Never give anything by mouth to an unconscious person. Give water to rinse the mouth.
- **Inhalation:** Move to fresh air.
- **Notes to physician:** Treat symptomatically.

5. Fire-fighting Measures

- **Suitable extinguishing media:** Any suitable media for extinguishing material supporting the fire
- **Special protective equipment for firefighters:** Wear self-contained breathing apparatus and protective suit.

6. Accidental Release Measures

- **Personal precautions:** Use personal protective equipment.
- **Methods for cleaning up:** Clean up all spills immediately, place in a suitable labelled container for waste disposal.

7. Handling and Storage

- **Handling:** Use good laboratory practices. Avoid inhaling vapors or mist.
Avoid contact with eyes, skin, or clothing. Wash hands following handling of material.
- **Storage :** Liquid Nitrogen storage. Store frozen cells at $\leq -130^{\circ}\text{C}$
- **Other precautions :** Explosion hazard; always follow biological safety precautions and liquid nitrogen safety rules.

8. Exposure Controls / Personal Protection

Exposure Limits

Engineering measures: Ensure adequate ventilation, especially in confined areas.

Personal Protective Equipment

Respiratory protection: In case of insufficient ventilation wear suitable respiratory equipment.

Hand protection: Impervious gloves.

Eye protection: Safety glasses with side-shields.

Skin and body protection: Lightweight protective clothing.

Hygiene measures: Handle in accordance with good industrial hygiene and safety practice.

Environmental exposure controls: Prevent product from entering drains.

9. Physical and Chemical Properties

General Information:

Form: Liquid

Appearance: Tinted Pink liquid in a cryovial

Odor: Odorless

Important Health Safety and Environmental Information:

Boiling point/range: No data available.

Melting poing/range: 0 °C.

Flash point: No data available.

Autoignition temperature: No data available.

Oxidising properties: No information available.

Water solubility: Soluble.

10. Stability and Reactivity

- **Stability:** Stable.
- **Materials to avoid:** Strong acids, acid chlorides, phosphorous halides, strong oxidizing agents, strong reducing agents.
- **Hazardous decomposition products:** Carbon monoxide, carbon dioxide, sulfur oxides.
- **Polymerization:** Hazardous polymerisation does not occur.

11. Toxicological Information

- **Acute toxicity**
- **Principle Routes of Exposure/Potential Health effects:**
 - Eyes: No information available
 - Skin: No information available
 - Inhalation: No information available
 - Ingestion: No information available
- **Specific effects**
 - Carcinogenic effects: No information available
 - Mutagenic effects: No information available
 - Reproductive toxicity: No information available
 - Sensitisation: No information available
- **Target Organ Effects:** No information available

12. Ecological Information

- **Ecotoxicity effects:** No information available
- **Mobility:** No information available
- **Biodegradation/ Bioaccumulation:** Inherently biodegradable/does not bioaccumulate

13. Disposal Considerations

Dispose of in accordance with local regulations.

14. Transport information

IATA:

- **Proper shipping name:** Not classified as dangerous in the meaning of transport regulations
- **Hazard Class:** No Information available.
- **Subsidiary Class:** No Information available.
- **Packaging Group:** No Information available.
- **UN-No:** No Information available.

15. Regulatory Information

- **Indication of danger :** Not hazardous.
- **R-phrase(s):** No Information available.
- **S-phrase(s):** No Information available.
- **International Inventories**

16. Other Information

The information provided herein was acquired by diligent search and/or investigation and the recommendations are based on prudent application of professional judgement. The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may present unknown hazards and should be used with caution. Since Novus Biologicals, LLC cannot control the actual methods, volumes, or conditions of use, the Company shall not be held liable for any damages or losses resulting from the handling or from contact with the product as described herein.

THE INFORMATION IN THIS MSDS DOES NOT CONSTITUTE A WARRANTY, EXPRESSED OR IMPLIED, INCLUDING ANY IMPLIED WARRANTY OF MERCHANTABILITY OR FITNESS FOR ANY PARTICULAR PURPOSE.

- **END OF SAFETY DATA SHEET**