

NutriFreez[®] D10 Cryopreservation Medium

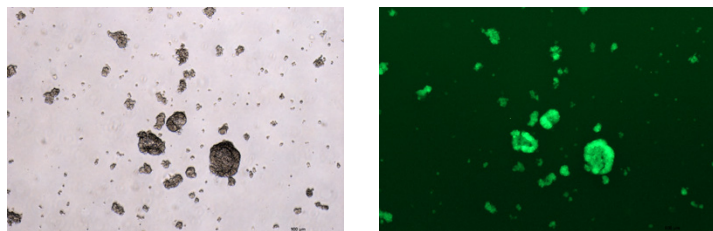
Powerful Cryopreservation
Media Optimized for Various
Cells and Tissues

NutriFreez[®] D10 Cryopreservation Medium is an optimized freezing solution designed and validated for the cryopreservation of various tissue and cell types, including but not limited to sensitive cell types such as hESCs, iPSCs, and MSCs. NutriFreez[®] D10 Medium maintains chemically defined and animal component-free conditions during cryopreservation, essential to maintaining consistency when culturing cells in a xeno-free system. NutriFreez[®] D10 Medium is ready-to-use and pre-formulated with DMSO, providing a protective environment for cells during the freezing, storage, and thawing process.



Cells preserved with NutriFreez[®] D10 Medium show excellent attachment (Figure 1) and maintain proper pluripotency marker expression after thawing, with superior results compared to both serum-containing freezing media, other serum-free solutions, and homebrew formulations¹.

Figure 1



Note. BGO1V/hOG cells (an Oct4-GFP reporter hES cell line) frozen in NutriFreez® D10 Medium and thawed into NutriStem® hPSC Medium on Matrigel. Images taken just 1 hour post-thaw show excellent survival and attachment of the hES cells, with high expression of Oct4 (green).

- High recovery post thaw
- Ready-to-use solution
- Serum-free and protein-free
- Chemically-defined
- cGMP-manufactured

Ordering Information

Cat. #	Product	Qty
05-713-1A	NutriFreez® D10 Cryopreservation Medium	500 mL
05-713-1B	NutriFreez® D10 Cryopreservation Medium	100 mL
05-713-1E	NutriFreez® D10 Cryopreservation Medium	50 mL
05-713-1C	NutriFreez® D10 Cryopreservation Medium	20 mL
05-713-1D	NutriFreez® D10 Cryopreservation Medium	10 mL
05-714-1A	NutriFreez® D10 Cryopreservation Medium, w/o phenol red	500 mL
05-714-1B	NutriFreez® D10 Cryopreservation Medium, w/o phenol red	100 mL
05-714-1E	NutriFreez® D10 Cryopreservation medium, w/o phenol red	50 mL
05-714-1D	NutriFreez® D10 Cryopreservation Medium, w/o phenol red	10 mL

Applicable Cell Types

- Human Embryonic Stem Cells
- Induced Pluripotent Stem Cells
- Human Mesenchymal Stem Cells
- Peripheral Blood Mononuclear Cells
- Human Endothelial Cells
- T cells, including Chimeric Antigen Receptor (CAR T) Cells and Tumor Infiltrating Lymphocytes (TILs)
- Neuron Cells
- Hybridomas
- CHO Cells
- Vero Cells
- Multiple mammalian cell lines: MRC-5, HEK293, HepG2, HeLa, BSC-1, BGM, 3T3, MA-10, BHK-21, B16-F10, MA-10

¹ Nishishita N, et al. An effective freezing | thawing method for human pluripotent stem cells cultured in chemically-defined and feeder-free conditions. *AJSC* 2015;4(1):38-49.

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

Israel

Biological Industries Israel Beit Haemek Ltd.
2511500 Kibbutz Beit Haemek
Phone: 972 4 9960595

 For more information, visit
www.sartorius.com