



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Thaw Medium 1
Catalog #: 60187-2
Lot #: 180323

Description

This liquid Cell Thawing Medium is optimized for thawing and plating select BPS Bioscience recombinant cell lines.

Size

500 ml. Sterile filtered (0.2 µm) and Mycoplasma-free

Quality

Each lot is tested for growth of recombinant cells. All lots are validated for low endotoxin levels (<1 EU/ml).

Storage and Stability

This medium is shipped on gel packs. Upon receipt, the medium should be stored in the dark at 4°C. If frozen during shipping, please move to 4°C for storage.. Protect from light. If the medium is warmed prior to use, do not exceed 37°C. When stored in the dark at 4°C, the product is stable at least 6 months from date of receipt.

Thawing Method

It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, transfer to a tube containing 10 ml of Thaw Medium 1, spin down cells, resuspend cells in pre-warmed Thaw Medium 1, transfer resuspended cells to T25 flask and culture at 37°C in a 7% CO2 incubator. At first passage switch to Thaw Medium 1 containing the correct amount of antibiotics (dependent upon the cell line). Cells should be split before they reach complete confluence. Reference individual cell line data sheets for more detailed thawing instructions.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Related Products

Product	Cat. #	Size
Thaw Medium 2	60184	100 ml
Thaw Medium 3	60186	100 ml
Thaw Medium 4	60181	100 ml
Thaw Medium 5	60182	100 ml
Thaw Medium 6	60183	100 ml
PD-1 Cell Growth Medium	60190	500 ml
PD-L1 Cell Growth Medium	60188	500 ml

Optimized For Use With

- CD155 (PVR) - HEK293 Recombinant Cell Line (Cat. # 60537)
- CD40/NF- κ B Reporter (Luc) - HEK293 Recombinant Stable Cell Line (Cat. # 60626)
- CD47 - HEK293 Cell Line (Cat. # 71249)
- CRE/CREB Reporter (Luc) – HEK293 Recombinant Cell Line (cAMP / PKA Signaling Pathway) (Cat. # 60515)
- GR-GAL4 Reporter (Luc)-HEK293 Recombinant Cell Line (Glucocorticoid Receptor Pathway) (Cat. # 60655)
- ISRE Reporter – HEK293 Recombinant Cell Line (JAK pathway) (Cat. # 60510)
- SLC5A5 – HEK293 Recombinant Cell line (Cat. #90333)
- SRE Reporter - HEK293 Recombinant Cell Line (ERK Pathway) (Cat. #60406)
- TDO-HEK293 Recombinant Cell line (Human) (Cat. #60534)
- TGF/SMAD Signaling Pathway SBE Reporter – HEK293 Cell Line (Cat. #60653)
- TRPC7-HEK293 Cell Line (Cat. #90030)
- NF- κ B reporter (Luc) - HEK293 Cell line (Cat. #60650)
- Notch Signaling Pathway Notch CSL Reporter – HEK293 Cell line (Cat. #60652)
- TCF/LEF reporter-HEK293 cell line (Wnt Signaling Pathway, Lithium-Dependent) (Cat. #60501)
- TRPM2-HEK293 Cell Line (Cat. #90331)
- hERG (Kv11.1) - HEK293 Recombinant Cell line (Cat. #60619)
- Empty vector control - HEK293 Recombinant Cell line (Cat. #90334)
- GAL4 Reporter (Luc)-HEK293 Cell Line (Cat. #60656)
- PDE7A-HEK293 Recombinant Cell line (Cat. #60407)
- PDE7B-HEK293 Recombinant Cell Line (Cat. #60412)
- hTMEM16A (ANO1) - HEK293 Recombinant Cell line (Cat. #90230)
- PDE7A/CRE Reporter - HEK293 Cell Line (Cat. #60413)
- hIDO1 - HEK293 Recombinant Cell Line (Cat. #60532)
- TMEM16B (ANO2) - HEK293 Recombinant Cell line (Cat. #90332)
- Hippo Pathway TEAD Reporter - MCF7 Recombinant Cell Line (Cat. #60618)
- IKCA1 (KCNN4) - HEK293 Recombinant Cell Line (Cat. #90330)
- Rat PDE7A - HEK293 Recombinant Cell line (Cat. #60408)
- hTRPC3-HEK293 Recombinant Cell line (Cat. #90130)
- PDE10A-HEK293 Recombinant Cell Line (Cat. #60410)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com