



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Thaw Medium 1**

### **Catalog #: 60187**

#### **Description**

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

#### **Size**

100 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. Do not freeze. 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 until the expiration date on the label (at least 2 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 5% CO<sub>2</sub> 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\*\*](mailto:info@bpsbioscience.com)

Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://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

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
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- $\kappa$ 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- $\kappa$ 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)

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

- 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\*\*](mailto:info@bpsbioscience.com)  
Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://www.bpsbioscience.com)