



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 3 Catalog #: 60186-2

Description

This liquid Cell Thaw Medium is optimized for thawing and plating BPS Bioscience's CHO cell-based recombinant cell lines, including

- TCR Activator - CHO Recombinant Cell line #60539
- TCR activator / PD-L1 - CHO Recombinant Cell line#60536
- D27 CHO-K1 Stable Recombinant Cell Line #60624
- NF- κ B Reporter (Luc) - CHO-K1 Recombinant Cell Line #60622

It is designed for use in a 37°C incubator with an atmosphere of 5% CO₂/95% air.

Note: once cells are well established, switch to an appropriate cell growth medium.

Size

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

Quality

Each lot is tested for growth of recombinant CHO 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 until 6 months from date of receipt).

Thawing Method

It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C waterbath, transfer to a tube containing 10 ml of Thaw Medium 3 **without antibiotics**, spin down the cells, and resuspend the cells in pre-warmed Thaw Medium 3 **without antibiotics**. Transfer resuspended cells to a T-25 flask and culture at 37°C in a CO₂ incubator overnight. The next day, replace the medium with Thaw Medium 3 **without antibiotics**, and continue growing culture in a CO₂ incubator at 37°C until the cells are ready to be split. Cells should be split before they reach complete confluence. At first passage, switch to Thaw Medium 3 **containing antibiotics**.

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

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
Thaw Medium 1	60187	100 ml
Thaw Medium 2	60184	100 ml
Thaw Medium 4	60181	100 ml
Thaw Medium 5	60182	100 ml
PD-1 Cell Growth Medium	60190	500 ml
PD-L1 Cell Growth Medium	60188	500 ml

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)