



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 2 Catalog #: 60184-1

DESCRIPTION:

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

- PD-1/NFAT Reporter-Jurkat cell #60535
- NFAT Reporter (Luc) – Jurkat Cell Line #60621
- NF-κB-Luciferase Reporter (Luc) - Jurkat Cell Line #60651
- ADCC Bioassay Effector Cell F variant (Low Affinity) #60540
- ADCC Bioassay Effector Cell V variant (High Affinity) #60541
- GITR / NF-κB-Luciferase Reporter (Luc) - Jurkat Cell Line #60546

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

SIZE:

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

QUALITY:

Each lot is tested for growth and functionality of luciferase reporter-Jurkat 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 water-bath, then transfer the entire contents of the vial to a tube containing 10 ml of Cell Thawing medium. Spin down the cells, remove supernatant and resuspend the cells in pre-warmed Cell Thawing Medium. Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO₂ incubator. Cells should be split before they reach 3x10⁶ cells/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

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
PD-1 Cell Growth Medium	#60189	500 ml
PD-L1 Cell Growth Medium	#60188	500 ml
PD-1/NFAT Reporter-Jurkat cell	#60535	2 vials
NFAT Reporter (Luc) – Jurkat Cell Line	#60621	2 vials
NF-κB Luciferase Reporter (Luc) - Jurkat Cell Line	#60651	2 vials
ADCC Bioassay Effector Cell F variant (Low Affinity)	#60540	2 vials
ADCC Bioassay Effector Cell V variant (High Affinity)	#60541	2 vials
Thaw Medium 1	#60187	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

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)