



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

Assay Medium 2A

Catalog #: 79621-2

DESCRIPTION:

This liquid Assay Medium is optimized for use in cellular assays in select BPS Bioscience's cell lines, including:

- ADCC Bioassay Effector Cell F variant (Low Affinity) – Jurkat Recombinant Cell Line #60540
- ADCC Bioassay Effector Cell V variant (High Affinity) – Jurkat Recombinant Cell Line #60541

SIZE:

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

FORMULATION:

RPMI, low IgG FBS 10%, 1% Penicillin/Streptomycin

QUALITY:

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.

RELATED PRODUCTS:

Product	Cat. #	Size
Thaw Medium 1	60187	100 ml
Thaw Medium 2	60184	100 ml
Growth Medium 2A	60190	500 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)