



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 2B

Catalog #: 79619-2

DESCRIPTION:

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

- Foxp3 Reporter (Luc) - Jurkat Recombinant Cell Line -#60628

SIZE:

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

FROMULATION:

Serum-free RPMI (Thermo Cat. #12045084), 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 2	60184	100 ml
Growth Medium 2B	79530	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

Please visit our website at: www.bpsbioscience.com