



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

## Growth Medium 2B Catalog #: 79530

### Description

This liquid Growth Medium is a complete medium optimized to culture select BPS Bioscience cell lines.

### Formulation

BPS Growth Media is formulated through BPS Thaw Media combined with antibiotics for maintenance of selective markers in the stable cell line.

- **Growth Medium 2B (BPS #79530):** Thaw Medium 2 (BPS #60184) plus 1 mg/ml of Geneticin (Thermo Fisher, #11811031).
  - **Thaw Medium 2 (BPS #60184):** RPMI 1640 medium (Thermo Fisher, #A1049101) supplemented with 10% FBS (Thermo Fisher, #26140079), 1% Penicillin/Streptomycin (Hyclone #SV30010.01).

### Quality

Sterile filtered (0.2  $\mu$ m) and Mycoplasma-free. Each lot is tested for growth of recombinant cells. All components are low endotoxin. FBS is ISIA-certified and sourced from the United States of America.

### Storage and Stability

Upon receipt, the medium should be stored in the dark at 4°C. Avoid repeated freeze/thaw cycles. Do not exceed 37°C. When stored in the dark at 4°C, the product is stable for six months after first thaw.

### Culture Protocol

For detailed thawing and cell culture protocol, see cell line data sheet.

### Optimized for Use With

Product	Cat. #	Size
IL-2 Jurkat Luciferase Reporter – Jurkat Cell Line	#60481	2 vials
NF-kB Luciferase Reporter – Jurkat Cell Line	#60651	2 vials
Foxp3 Reporter (Luc) – Jurkat Recombinant Cell Line	#60628	2 vials

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)