



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Growth Medium 2A**

### **Catalog #: 60190**

#### **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 2A (BPS #60190):** Thaw Medium 2 (BPS #60184) plus 1 mg/ml of Geneticin (Thermo Fisher, #11811031), and 200 µg/ml of Hygromycin B (Thermo Fisher, #10687010).
  - **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 µm) 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.

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Optimized for Use With**

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
PD-1 / NFAT Reporter - Jurkat Recombinant Cell Line	#60535	2 vials
LAG3 / NFAT Reporter - Jurkat Recombinant Cell Line	#71278	2 vials
ADCC Bioassay Effector Cell V variant (High Affinity) Jurkat Recombinant Cell Line	#60541	2 vials
ADCC Bioassay Effector Cell F variant (Low Affinity) Jurkat Recombinant Cell Line	#60540	2 vials
NFAT Reporter (Luc) – Jurkat Recombinant Cell Line	#60621	2 vials
HVEM/NF- $\kappa$ B Reporter Jurkat Recombinant Cell Line	#79310	2 vials
TIGIT / NFAT Reporter - Jurkat Cell Line	#60538	2 vials
ADCP Bioassay Effector Cell Fc $\gamma$ RIIa (H variant) NFAT Reporter-Jurkat	#71273	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)