



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 8A**

### **Catalog #: 79653**

#### **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 8A (BPS #79653):** Thaw Medium 8 (BPS #79652) plus 1 µg/ml of Puromycin (Takara, #631306).
  - **Thaw Medium 8 (BPS #79652):** RPMI 1640 medium (Thermo Fisher, #A1049101) supplemented with 10% heat-inactivated FBS (Thermo Fisher, #10082147), 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.

#### **Optimized for Use With:**

- NF- κB Reporter (Luc) – THP-1 Cell Line - #79645

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)