



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Growth Medium 10A Catalog #: 79835**

### **Description**

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

### **Formulation**

BPS Growth Medium 10A is formulated through BPS Growth Medium combined with antibiotics for maintenance of selective markers in the stable cell line.

- **Growth Medium 10A (BPS Bioscience, #79835):** Growth medium with 400 µg/ml of Geneticin (Life technologies, #11811031) to ensure recombinant expression.
  - **Thaw Medium 10 (BPS Bioscience, #79704):** RPMI 1640 medium supplemented with 10% FBS, 1 MM Sodium pyruvate, 1 % Non-essential amino acids, 1 % Penicillin/streptomycin.

### **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:**

- Firefly Luciferase – RPMI 8226 Recombinant Cell Line - #79834

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** 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)