



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)

## **BPS Medium 7 Catalog #: 79260**

### **Description**

This liquid BPS Medium is optimized for use with select BPS reporter kits.

### **Size**

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

### **Quality**

Each lot is tested for growth of recombinant cells. 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 for 6 months from date of receipt.

### **Growth Method**

See individual protocols for detailed instructions.

### **Related Products**

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
Thaw Medium 1	60187	100 ml
Thaw Medium 2	60184	100 ml
Thaw Medium 3	60186	100 ml
Thaw Medium 4	60181	100 ml
Thaw Medium 5	60182	100 ml
Thaw Medium 6	60183	100 ml

### **Optimized For Use With**

- Myc Reporter Kit (BPS Cat. #60519)
- Transfection Collection™ : Myc Transient Pack (Myc Signaling Pathway) (BPS Cat. #79284)

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)