



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)

## **Thaw Medium 7**

### **Catalog #: 60185**

#### **Description**

This liquid Cell Thawing Medium is optimized for thawing and plating select BPS Bioscience recombinant cell lines, including:

- Myc Reporter (Luc)-HCT116 Recombinant Cell Line (Myc Signaling Pathway) (BPS Bioscience #60520)
- NF- $\kappa$ B Reporter (Luc) -HCT116 Recombinant Cell Line (BPS Bioscience #60623)

#### **Formulation**

McCoy's 5A medium (Hyclone #SH30200.01) with 10% FBS (Thermo Fisher, Cat. #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.

#### **Thawing Protocol**

For detailed thawing 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 **1.858.202.1401** 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)