



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)

## **Assay Medium 1A**

### **Catalog #: 79805**

#### **DESCRIPTION:**

This liquid Assay Medium is optimized for use in cellular assays in select BPS Bioscience's cell lines, including:

- TrkA / SRE Reporter – HEK293 Recombinant Cell Line (#79798)

#### **FORMULATION:**

Opti-MEM (Thermo Fisher, #31985070) plus 1% Penicillin/Streptomycin (Hyclone, #SV30010.01).

#### **QUALITY:**

All lots are validated for low endotoxin levels (<1 EU/ml).

#### **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.

#### **RELATED PRODUCTS:**

<b>Product</b>	<b>Cat. #</b>	<b>Size</b>
TrkA / SRE Reporter – HEK293 Recombinant Cell Line	79798	2 vials

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

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