



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

## Growth Medium 3E Catalog #: 79553

### 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 3E (BPS #79553):** Thaw Medium 3 (BPS #60186) plus 100 µg/ml of Hygromycin B (Thermo Fisher, #10687010).
  - **Thaw Medium 3 (BPS #60186):** Ham's F-12 medium (Hyclone # SH30526.01) supplemented with 10% FBS (Thermo Fisher, #26140079), 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

Product	Cat. #	Size
CD27 CHO-K1 Stable Recombinant Cell Line	#60624	2 vials
HVEM - CHO Recombinant Cell Line	#79297	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)