



6405 Mira Mesa Blvd Ste 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## Growth Medium 3D Catalog #: 79539

### 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 3D (BPS #79539):** Thaw Medium 3 (BPS #60186) plus 1 mg/ml Geneticin (Thermo Fisher, #11811031).
  - **Thaw Medium 3 (BPS #60186):** F-12K Medium (Kaighn's Modification of Ham's F-12 Medium) supplemented with 10% FBS (Thermo Fisher, #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.

### Culture Protocol

For detailed thawing and cell culture protocol, see cell line data sheet.

### Optimized for Use With

Product	Cat. #	Size
LIGHT-CHO Recombinant Cell Line	#79262	2 vials
PD-L1 - CHO Recombinant Cell Line	#60543	2 vials
NF- $\kappa$ B Reporter (Luc) - CHO-K1 Recombinant Cell Line	#60622	2 vials
CD137L (4-1BBL) CHO-K1 Recombinant Cell Line	#60523	2 vials
GITRL CHO-K1 Recombinant Cell Line	#60547	2 vials
CD27 CHO-K1 Stable Recombinant Cell Line	#60624	2 vials
Transfection Collection™ : NF- $\kappa$ B Reporter Cellular Assay Pack (CHOK1)	#79325	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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)