



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Growth Medium 3G Catalog #: 79882

Description

This liquid Growth Medium is a complete medium optimized to culture select BPS Bioscience cell lines and lentivirus.

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 3G (BPS Bioscience #79882):** Thaw Medium 3 (BPS Bioscience #60186) plus 1000 µg/ml Geneticin (Thermo Fisher, #11811031) and 5 µg/ml Puromycin (InvivoGen, #ant-pr-1).
 - **Thaw Medium 3 (BPS Bioscience #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 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
FcGR11A (CD16a) Lentivirus	#79876	2 vials
FcER1G Lentivirus	#79878	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

Please visit our website at: www.bpsbioscience.com