



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Growth Medium 5B Catalog #: 79541

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 5B (BPS #79541):** Thaw Medium 5 (BPS #60182) plus 500 µg/ml of Geneticin (Thermo Fisher, #11811031).
 - **Thaw Medium 5 (BPS #60182):** DMEM medium (Hyclone #SH30243.01) supplemented with 10% Calf Serum (CS) (Hyclone # SH30072.03) (not FBS), 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. Calf Serum 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

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
Gli Reporter – NIH3T3 Recombinant Cell line (Hedgehog Pathway)	#60409	2 vials

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com