



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 4B

Catalog #: 79554

Description

This liquid Growth Medium is a complete medium optimized to culture select BPS Bioscience cell lines. Follow the culturing instructions on the designated protocol.

Size

500 ml. Sterile filtered (0.2 µm) and Mycoplasma-free

Formulation

DMEM/F12 (1:1) (Hyclone # SH30271.01) supplemented with 10% FBS (Life Technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone # SV30010), 10 µg/ml Blasticidin (Life Technologies # R210-01), 200 ug /ml Zeocin (Invivogen # ant-zn-1p), and 100 ug/ml Hygromycin B (Hyclone #SV30070.01)

FBS is ISIA-certified and sourced from BSE-free regions.

Quality

Each lot is tested for growth of a select cell line. All lots are validated for low endotoxin levels (<1 EU/ml).

Storage and Stability

This medium is shipped on gel packs. Upon receipt, the medium should be stored in the dark at 4°C. If frozen during shipping, please move to 4°C for storage. Protect from light. If the medium is warmed prior to use, do not exceed 37°C. When stored in the dark at 4°C, the product is stable for 6 months from date of receipt.

Optimized for Use With

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
Nav1.8/β2 - HEK293 Recombinant Cell Line	#60521	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