



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 2F Catalog #: 79669

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 2F (BPS #79669):** Thaw Medium 2 (BPS #60184) plus 1 mg/ml Geneticin (Thermo Fisher, #11811031), 0.5 µg/ml Puromycin Dihydrochloride (Thermo Fisher, #A1113803).
 - **Thaw Medium 2 (BPS #60184):** RPMI 1640 medium (Thermo Fisher, #A1049101) 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

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
ICOS NFAT - Jurkat Cell Line	#79668	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

Please visit our website at: www.bpsbioscience.com