



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Growth Medium 1F**

### **Catalog #: 79540**

#### **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.

#### **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 1F (BPS #79540):** Thaw Medium 1 (BPS #60187) plus 100 µg/ml Hygromycin B (Thermo Fisher, #10687010).
  - **Thaw Medium 1 (BPS #60187):** MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Invitrogen #26140-079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 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
CD155 (PVR) - HEK293 Recombinant Cell Line	#60537	2 vials
PD-1 - HEK293 Recombinant Cell Line	#60680	2 vials
CRE/CREB Reporter (Luc) – HEK293 Recombinant Cell Line (cAMP / PKA Signaling Pathway)	#60515	2 vials
CTLA4 - HEK293 Recombinant Cell Line	#60681	2 vials
GITR - HEK293 Recombinant Cell Line	#79092	2 vials
TMEM16B (ANO2) - HEK293 Recombinant Cell line	#90332	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](mailto:info@bpsbioscience.com)

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