



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

BPS Medium 8 Catalog #: 79385

Description

This liquid BPS Medium is optimized for use with the CSF1R/SRE Reporter Kit (BPS Bioscience, #79379).

Size

100 ml. Sterile filtered (0.2 μ m) and Mycoplasma-free

Quality

Each lot is tested for growth of recombinant cells. 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 the date of receipt.

Growth Method

See individual protocols for detailed instructions.

Related Products

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
Thaw Medium 1	60187	100 ml
Thaw Medium 2	60184	100 ml
Thaw Medium 3	60186	100 ml
Thaw Medium 4	60181	100 ml
Thaw Medium 5	60182	100 ml
Thaw Medium 6	60183	100 ml
PD-1 Cell Growth Medium	60190	500 ml
PD-L1 Cell Growth Medium	60188	500 ml

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)



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

Optimized For Use With

- AP1 Reporter Kit (JNK Pathway) (Cat. #60612)
- ARE Reporter Kit (Nrf2 Antioxidant Pathway) (Cat. #60514)
- CRE/CREB Reporter Kit (cAMP/PKA Signaling Pathway) (Cat. #60611)
- NF- κ B Reporter Kit (NF- κ B Signaling Pathway) (Cat. #60614)
- Notch1/CSL Reporter Kit (Notch Signaling Pathway) (Cat. #60509)
- SBE Reporter Kit (TGF β /SMAD signaling pathway) (Cat. #60654)
- SRE Reporter Kit (MAPK/ERK Signaling Pathway) (Cat. #60511)
- TCF/LEF Reporter Kit (Wnt Signaling Pathway) (Cat. #60500)

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)