



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 1 Catalog #: 79259

Description

This liquid BPS Medium is optimized for use with many BPS Reporter Kits.

Size

100 ml. Sterile filtered (0.2 µm) 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

Optimized For Use With

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

- 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)