



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

Dear customer:

Thank you for choosing BPS Bioscience for your assay and screening needs! Here are some helpful tips when shipping us compounds.

1. Compound Preparation:

- i. (preferred) Prepare at least 50 μ l per compound, at a concentration of 10 mM in DMSO
- ii. Powder preparations are acceptable as long as the MW of the compounds are included to ensure accurate dilutions are made.
- iii. If sending a powder, please send 1-3 mg, depending on the MW of the compound.

2. Compound Submission Form:

- i. You can find this online at <http://bpsbioscience.com/compound-submissions>.
- ii. *For all:* fill in column B and check either column C (for IC/EC50) or column D (inhibition screen)
- iii. *Assay information:* For IC50/inhibition screens fill in columns E-I. For cell-based assays fill in columns J-N
- iv. *Compound information:* For liquid stocks fill in columns O-R. For powder stocks fill in columns S-W
- v. *Additional notes* can be placed in column X

3. Shipping:

- i. Please ship to the following address:

BPS Bioscience, Inc.
ATTN: Andrew Newman
6042 Cornerstone Ct. W., Suite B
San Diego, CA 92121 USA
Tel #: (858) 829-3082

- ii. Please include the Estimate # or printed Compound submission form with your shipment for optimal processing
- iii. For international shipments: Please also include origin, destination, and package contents in English. Many couriers require a "Declaration of Biological Shipment." Please follow the instructions and fill this out appropriately.
- iv. After shipping, please email service-orders@bpsbioscience.com with your tracking number so we can monitor your shipment.

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