



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### ***SHP-2 Substrate (DiFMUP)***

**Catalog #: 79769**

**Description:** The fluorogenic SHP-2 substrate, DiFMUP, is added to the mixture of SHP-2 (BPS Bioscience #79018) and SHP-2 Activating Peptide (BPS Bioscience #79319). DiFMU fluorophore is released and can then be measured using a fluorescence reader.

**Background:** DiFMUP is a fluorinated 4-methylumbelliferyl phosphate (MUP) that is ideal as a substrate for fast, sensitive phosphatase assays, even at low pH. DiFMUP has an excitation/emission maxima of ~358/450 nm.

**Concentration:** 1.0 mM

**Supplied as:** DMSO solution

**Size:** 50 µl

**Storage:** -80°C

**Stability:** Stable for up to 6 months at -80°C when stored as directed. Avoid freeze/thaw cycles.

**Reference:** Welte, S., *et al. Anal Biochem.* 2005 Mar 1;**338(1)**:32-8.

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](mailto:info@bpsbioscience.com)

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