



6405 Mira Mesa Blvd Ste. 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

**Data Sheet**  
**Fluorogenic SHP1 (PTPN6) Assay Kit**  
Catalog #79831  
Size: 96 reactions

**BACKGROUND:** SHP1 (also known as PTPN6) is preferentially expressed in a variety of hematopoietic cells, and is an early response gene in lymphokine-stimulated cells. The noncatalytic N-terminus of this phosphatase can interact with MAP kinases and negatively regulates ERK2 and p38 MAP-kinases activity. The PTPN6 is involved in the regulation of T cell antigen receptor (TCR) signaling, which is thought to function through dephosphorylating molecules related to the MAP kinase pathway.

**DESCRIPTION:** The *Fluorogenic SHP1 Assay Kit* is designed to measure SHP1 activity for screening and profiling applications, in a homogeneous assay with no time-consuming washing steps. The SHP1 assay kit comes in a convenient 96-well format, with purified SHP1 enzyme, fluorogenic substrate, and PTP assay buffer for 100 enzyme reactions. This kit is developed to identify inhibitors of the regulatory region of SHP1.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
100342	Recombinant Human SHP1 (221-523)	≥1 µg	-80°C	<b>Avoid freeze/ thaw cycles!</b>
	0.5 mM PTP Substrate (DiFMUP)	50 µl	-80°C	
79716	5X PTP Assay Buffer	3 ml	-20°C	
	0.5 M DTT	200 µl	-20°C	
79685	Black, low binding black microtiter plate	1	Room Temperature	

**MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

Fluorescent microplate reader capable of reading exc/em=360nm/460nm

**APPLICATIONS:** Great for studying enzyme kinetics and HTS applications.

**STABILITY:** One year from date of receipt when stored as directed.

**REFERENCE(S):**

Gee, K.R., *et al.*, *Anal Biochem*, 1999 Aug 15; **273(1)**:41-8.

Kilgore, N.E., *et al.*, *J. Immunol.* 2003, **170**:4891-4895.

Duchesne, C., *et al.* *J. Biol. Chem* 2003, **278**:14274-14283.

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd Ste. 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## **ASSAY PROTOCOL:**

***All samples and controls should be tested in duplicate.***

- 1) Prepare 1X assay buffer with 1 mM DTT from 5X assay buffer. For example, add 200  $\mu$ l 5X assay buffer and 2  $\mu$ l 0.5 M DTT to 798  $\mu$ l distilled H<sub>2</sub>O to make 1 ml 1X assay buffer.
- 2) Thaw SHP1 on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot SHP1 into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. *Note: SHP1 enzyme is sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 3) Dilute SHP1 in 1x assay buffer at 0.15 ng/ $\mu$ l (3 ng per reaction).
- 4) Add 20  $\mu$ l diluted SHP1 enzyme solution to wells designated "Positive Control" and "Test Sample." Add 20  $\mu$ l 1X assay buffer to "Blank" wells.
- 5) Prepare the inhibitor solution.

If the inhibitor compound is dissolved in DMSO, make a 100-fold higher concentration of the compound than the highest concentration you want to test in DMSO. Then make a 10 fold dilution in water (at this step the compound concentration is 10-fold higher than the final concentration in 10% DMSO). If you want to an run IC<sub>50</sub> or test lower concentrations of the compound, prepare a series of further dilutions in water containing 10% DMSO (final concentration of the DMSO will be 1% in all samples).

If the inhibitor compound is dissolved in water, make a solution of the compound 10-fold higher than the final concentration in water.

- 6) Add 5  $\mu$ l inhibitor solution to each well designated "Test Sample." Add 5  $\mu$ l of 10% DMSO in water (inhibitor buffer) to "Blank" and "Positive Control" wells.
- 7) Incubate the enzyme with the compound solution at room temperature for 30 minutes.

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd Ste. 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

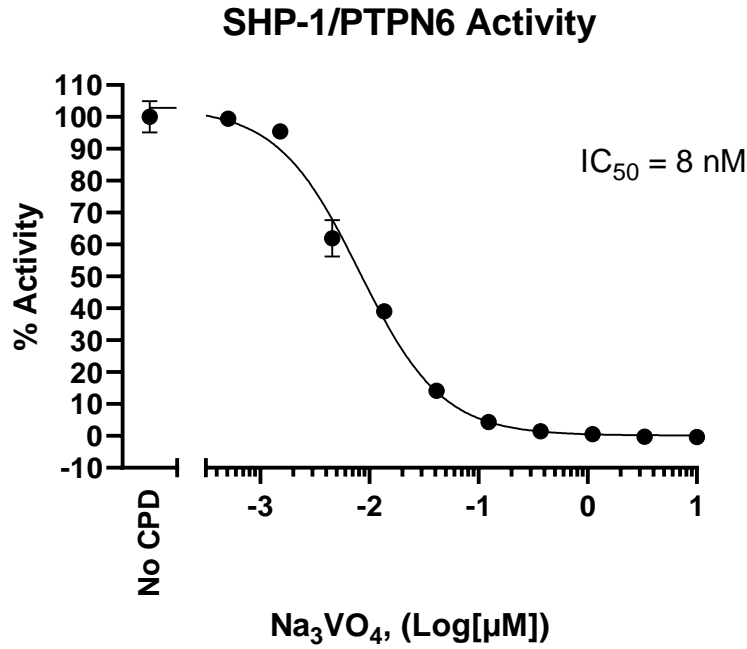
Component	Positive Control	Test Sample	Blank
SHP1 (0.15 ng/μl)	20 μl	20 μl	20 μl
Test Inhibitor	-	5 μl	-
10% DMSO in Water (inhibitor buffer)	5 μl	-	5 μl
Incubate at room temperature for 30 minutes			
Substrate solution	25 μl	25 μl	25 μl
<b>Total</b>	<b>50 μl</b>	<b>50 μl</b>	<b>50 μl</b>

- 8) Prepare the substrate solution: N wells × (24.5 μl 1X assay buffer (with DTT) + 0.5 μl 0.5 mM PTP Substrate).
- 9) Add 25 μl of the substrate solution to each well (Final concentration of the PTP substrate in a 50 μl reaction is 5 μM).
- 10) Incubate at room temperature for 30 minutes. Measure the fluorescence intensity in a microtiter plate-reading fluorimeter capable of excitation at a wavelength 360 nm and detection of emission at a wavelength 460 nm. The fluorescence intensity can also be measured kinetically. "Blank" value is subtracted from all other values.

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

**Example of Assay Results:**



SHP1 enzyme activity, measured using the *Fluorogenic SHP1 Assay Kit* (BPS Bioscience #79831). Fluorescence intensity was measured using a Tecan fluorescent microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

**Related Products**

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
SHP-1, His-Tag	100342	20 $\mu$ g
SHP-1 (PTPN6)	30021	20 $\mu$ g
5X PTP Assay Buffer	79716	10 ml

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**  
 Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)