

**Data Sheet**  
***SHP-2 (catalytic domain) Homogeneous Assay Kit***  
**Catalog #79317**  
**Size: 96 reactions**

**DESCRIPTION:** Mammalian PTPases can be subdivided into 1 of 2 broad categories: transmembrane receptor PTPases and intracellular PTPases. SHP-2 (PTPN11) is one of the 2 closely related mammalian intracellular PTPases whose sequences encode 2 tandem SRC homology 2 (SH2) domains that are located at the amino-terminal side of a single PTPase catalytic domain. This PTP is widely expressed in most tissues and plays a regulatory role in various cell signaling events that are important for a diversity of cell functions, such as mitogenic activation, metabolic control, transcription regulation, and cell migration.

The SHP-2 (catalytic domain) Homogeneous Assay Kit is a complete assay system designed to measure SHP-2 activity for screening and profiling applications. It comes in a convenient 96-well format, with all the reagents necessary for 100 fluorescent SHP-2 activity measurements. In addition, the kit includes purified SHP-2 enzyme and SHP-2 inhibitor, orthovanadate, for use as a positive and negative control. Using this kit, only one simple step on a microtiter plate is needed to analyze the SHP-2 activity level. The fluorogenic substrate, DiFMUP, is incubated with purified SHP-2 and the enzymatic activity releases DiFMU fluorophore that can then be measured using a fluorescence reader.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
30022	Recombinant Human SHP-2	1 µg	-80°C	<b><i>Avoid freeze/ thaw cycles!</i></b>
79769	1mM SHP-2 Substrate (DiFMUP)	50 µl	-80°C	
79626	5X SHP-2 Assay Buffer	3 ml	-20°C	
	1mM Na <sub>3</sub> VO <sub>4</sub>	20 µl	-80°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):**

Chai, J., et al., Cell, 2001 Mar 9;104(5):769-80.

Denault, JB., and Salvesen, GS., J. Biol. Chem., 2003 Sep 5;278(36):34042-50.

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)

**ASSAY PROTOCOL:**

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

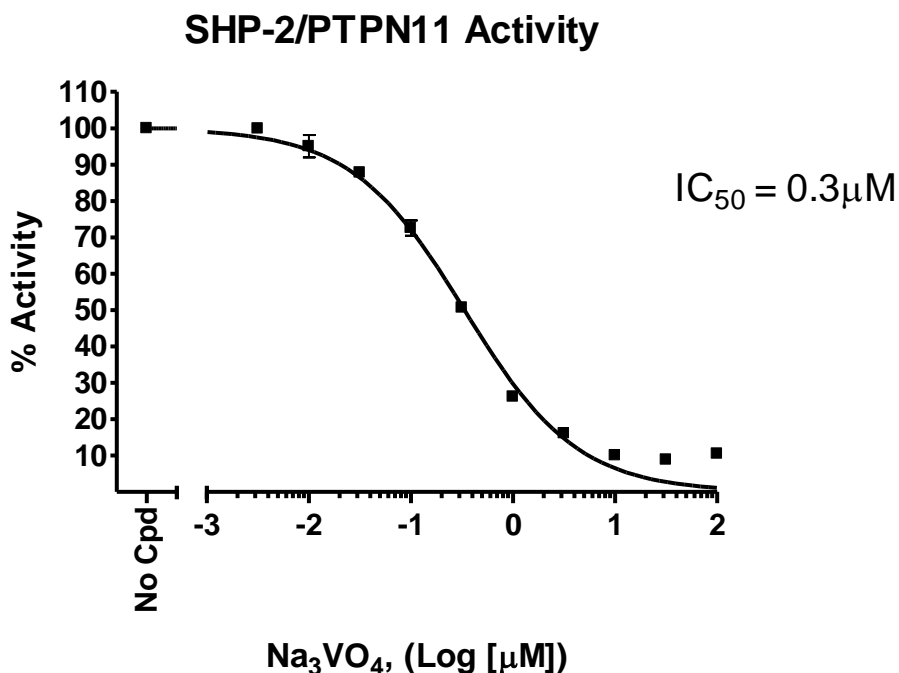
- 1) Prepare 1X assay buffer from 5X assay buffer. For example, add 200 $\mu$ l 5X assay buffer to 800 $\mu$ l H<sub>2</sub>O to make 1ml 1X assay buffer.
- 2) Prepare the master mixture: N wells  $\times$  (19.5  $\mu$ l H<sub>2</sub>O, 5  $\mu$ l 5X SHP-2 assay buffer + 0.5  $\mu$ l 1 mM SHP-2 substrate (DiFMUP)).
- 3) Add 25  $\mu$ l of master mixture to each well (The final substrate concentration in a 50  $\mu$ l reaction is 10  $\mu$ M).
- 4) Prepare the inhibitor solution that is 10-fold higher than the final concentration.
- 5) Add 5  $\mu$ l of the inhibitor solution to the well designed with "Test Sample". Add 5  $\mu$ l of the inhibitor buffer (without inhibitor) to the wells designed with "Blank" and "Positive Control". Add 5  $\mu$ l of control compound, Na<sub>3</sub>VO<sub>4</sub>, to the well designed with "Negative Control".
- 6) Thaw SHP-2 on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot SHP-2 into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. *Note: SHP-2 enzyme is sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 7) Dilute SHP-2 in 1x assay buffer at 0.2 ng/ $\mu$ l (4 ng per reaction).
- 8) Add 20  $\mu$ l diluted SHP-2 solution to the wells designed with "Positive Control", "Test Sample" and Negative Control".
- 9) 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. "Blank" value is subtracted from all other values. You can also measure the fluorescence intensity kinetically.

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)

Component	Positive Control	Test Sample	Negative Control	Blank
H <sub>2</sub> O	19.5 µl	19.5 µl	19.5 µl	40 µl
5X assay buffer	5 µl	5 µl	5 µl	5 µl
SHP-2 Substrate (DiFMUP)	0.5 µl	0.5 µl	0.5 µl	-
Test Inhibitor	-	5 µl	-	-
Control Inhibitor	-	-	5 µl	-
Inhibitor Buffer (no inhibitor)	5 µl	-	-	5 µl
SHP-2 (0.2 ng/µl)	20 µl	20 µl	20 µl	-
<b>Total</b>	<b>50 µl</b>	<b>50 µl</b>	<b>50 µl</b>	

**Example of Assay Results:**



SHP-2 enzyme activity, measured using the *SHP-2 (catalytic domain) Homogeneous Assay Kit*, BPS Bioscience #79317. 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)

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)