



6042 Cornerstone Court W, Ste 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**  
**SPHK1 Assay Kit**  
Catalog # 78026  
96 Reactions

**BACKGROUND:** SPHK1 is a lipid enzyme that catalyzes the phosphorylation of sphingosine to form sphingosine 1-phosphate (SPP). It has been implicated in tumor progression through cell proliferation and motility. High levels of SPHK1 have been associated with increased mortality in various forms of human cancer.

**DESCRIPTION:** The *SPHK1 Assay Kit* is designed to measure SPHK1 activity for screening and profiling applications using Kinase-Glo<sup>®</sup> MAX as a detection reagent. The *SPHK1 Assay Kit* comes in a convenient 96-well format, with enough purified SPHK1, Sphingosine, ATP, and kinase assay buffer for 96 enzyme reactions.

**COMPONENTS:**

Catalog #	Reagent	Amount	Storage	
40610	SPHK1, His-tag*	>1 µg	-80°C	<b>Avoid multiple freeze/ thaw cycles!</b>
79334	5x Kinase assay buffer	1.5 ml	-20°C	
79686	ATP (500 µM)	100 µl	-20°C	
	Sphingosine (1 mM)	100 µl	-20°C	
79696	96-well plate, white	1	RT	

\*Excess material has been provided for ease of retrieval from the vial.

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

Kinase-Glo MAX (Promega #V6071)  
Dithiothreitol (DTT, 0.5 M)  
Microplate reader capable of reading luminescence  
Adjustable micropipettor and sterile tips  
30°C incubator

**APPLICATIONS:** Useful for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

**STABILITY:** Up to 6 months when stored as recommended.

**REFERENCES:**

1. Shida, D., *et al.* 2008. "Targeting SphK1 as a new strategy against cancer." *Current Drug Targets* **9(8)**: 662-673.
2. Xia, J., *et al.* 2012. "miR-124 inhibits cell proliferation in gastric cancer through down-regulation of SPHK1." *The Journal of Pathology* **227(4)**: 470-480.

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)

**ASSAY PROTOCOL:**

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

- 1) Thaw **5x Kinase assay buffer**, **ATP (500  $\mu\text{M}$ )**, and **Sphingosine (1 mM)**. Add 30  $\mu\text{l}$  of 0.5 M DTT to **5x Kinase assay buffer**.
- 2) Prepare the master mixture (25  $\mu\text{l}$  per well): N wells x (5  $\mu\text{l}$  **5x Kinase assay buffer** + 1  $\mu\text{l}$  **ATP (500  $\mu\text{M}$ )** + 1  $\mu\text{l}$  **Sphingosine (1 mM)** + 18  $\mu\text{l}$  distilled water). Add 25  $\mu\text{l}$  to every well.

	Positive Control	Test Inhibitor	Blank
5x Kinase assay buffer	5 $\mu\text{l}$	5 $\mu\text{l}$	5 $\mu\text{l}$
ATP (500 $\mu\text{M}$ )	1 $\mu\text{l}$	1 $\mu\text{l}$	1 $\mu\text{l}$
Sphingosine (1 mM)	1 $\mu\text{l}$	1 $\mu\text{l}$	1 $\mu\text{l}$
Water	18 $\mu\text{l}$	18 $\mu\text{l}$	18 $\mu\text{l}$
Test Inhibitor	-	5 $\mu\text{l}$	-
Inhibitor buffer (10% DMSO in water)	5 $\mu\text{l}$	-	5 $\mu\text{l}$
1x Kinase buffer	-	-	20 $\mu\text{l}$
SPHK1, His-tag (0.19 ng/ $\mu\text{l}$ )	20 $\mu\text{l}$	20 $\mu\text{l}$	-
<b>Total</b>	<b>50 <math>\mu\text{l}</math></b>	<b>50 <math>\mu\text{l}</math></b>	<b>50 <math>\mu\text{l}</math></b>

- 3) Add 5  $\mu\text{l}$  of Inhibitor solution of each well labeled as "Test Inhibitor." For the "Positive Control" and "Blank," add 5  $\mu\text{l}$  of the same solution without inhibitor (Inhibitor buffer, usually 10% DMSO in water). *Note: Final DMSO concentration must be  $\leq 1\%$ . Higher DMSO levels can significantly decrease the enzyme activity. For example, to test an inhibitor at 10  $\mu\text{M}$  that is dissolved in 100% DMSO, dilute 1 mM inhibitor with water to make a 100  $\mu\text{M}$  inhibitor in 10% DMSO(aq). Then, add 5  $\mu\text{l}$  of the 100  $\mu\text{M}$  solution into the 50  $\mu\text{l}$  assay to make a 1% DMSO concentration in the final reaction mixture.*
- 4) Prepare 3 ml of **1x Kinase assay buffer** by mixing 600  $\mu\text{l}$  of **5x Kinase assay buffer** with 2400  $\mu\text{l}$  water. 3 ml of **1x Kinase assay buffer** is sufficient for 100 reactions.
- 5) To the wells designated as "Blank," add 20  $\mu\text{l}$  of **1x Kinase assay buffer**.
- 6) Thaw **SPHK1, His-tag** on ice. Upon first thaw, briefly spin tube containing material to recover full content of the tube. Calculate the amount SPHK1, His-tag required for the assay and dilute enzyme to 0.19 ng/ $\mu\text{l}$  with **1x Kinase assay buffer**. Store remaining undiluted material in aliquots at  $-80^{\circ}\text{C}$ . *Note: SPHK1,*

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**

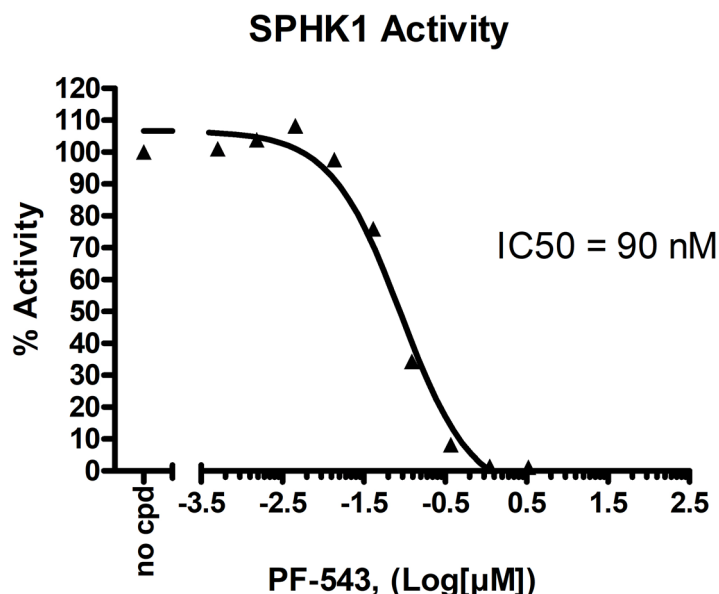
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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

*His-tag is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted material.*

- 7) Initiate reaction by adding 20  $\mu$ l of diluted **SPHK1**, **His-tag** to the wells designated "Positive Control" and "Test Inhibitor Control." Incubate at 30°C for 45 minutes.
- 8) Thaw Kinase-Glo Max reagent.
- 9) After the 45-minute reaction, add 50  $\mu$ l of Kinase-Glo Max reagent to each well. Cover plate with aluminum foil and incubate the plate at room temperature for 15 minutes.
- 10) Measure luminescence using the microplate reader. The value of the "Blank" reading should be subtracted from all other measurements.

#### Example of Assay Results:



Inhibition of SPHK1, His-tag by PF-543, measured using the SPHK1 assay kit (BPS Bioscience #78026). *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@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: [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:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<b><u>Product Name</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
Sphingosine kinase 1, His-tag	40610	20 µg
5x Kinase assay buffer	79334	10 ml
ATP (500 µM)	79686	200 µl
Protein Tyrosine Kinase Substrate (poly-Glu,Tyr 4:1)	40217	1 mg
Sphingosine kinase 2, His-tag	40611	20 µg
Sphingosine kinase 1, His-tag	40610	20 µg
Sphingosine kinase 2 (long), His-tag	40612	10 µg
Mouse Sphingosine kinase 1a, His-tag	40613	10 µg
Mouse Sphingosine kinase 2, His-tag	40614	10 µg

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)