

## Data Sheet

### SPHK2

Human, recombinant, N-terminal HIS tag  
**Catalog #:** 40611  
**Lot#:** 100219      **Conc.:** 0.5 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
124 mM NaCl, 2.4 mM KCl, 225 mM  
imidazole, 10% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80 °C

**References:**

1. Zemann, B. *et al.*, *Blood*. **107(4)**:1454-8 (2006)
2. Maceyka, M. *et al.*, *J Biol Chem*. **280(44)**:37118-29 (2005)

**Description:** Human Sphingosine kinase 2 (GenBank Accession No. AF245447), full length with N-terminal His tag, MW=69.5 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 20 U/μg. Sphingosine kinase activity was determined using D-erythro-sphingosine and ATP as substrates. A unit of sphingosine kinase activity is defined as the amount of enzyme required to produce 1 pmol of S1P/min at 37°C. Assay condition: 50 mM HEPES, pH 7.4, 150 mM NaCl, 5 mM MgCl<sub>2</sub>, 1 mM DTT, 3 μM Na-orthovanadate, 0.5 mM ATP, 4 μM D-erythro-sphingosine, and 0.75 μg/ml sphingosine kinase 2 at 37°C for 30 min. ATP reduction is detected using Kinase-Glo<sup>®</sup> Luminescent Kinase Assay Platform (Promega Corporation, Madison, WI).

**Application:**

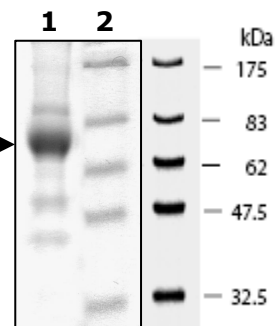
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
15 μg SPHK2  
**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 69.5 kDa.  
**Purity:** ≥80%



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](mailto:info@bpsbioscience.com)

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