

Data Sheet

SPHK2, His-tag (Mouse)

Mouse recombinant, N-terminal His-tag

Catalog #: 40614

Lot #: 190523-2

Conc.: 0.57 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 3mM DTT and 50% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

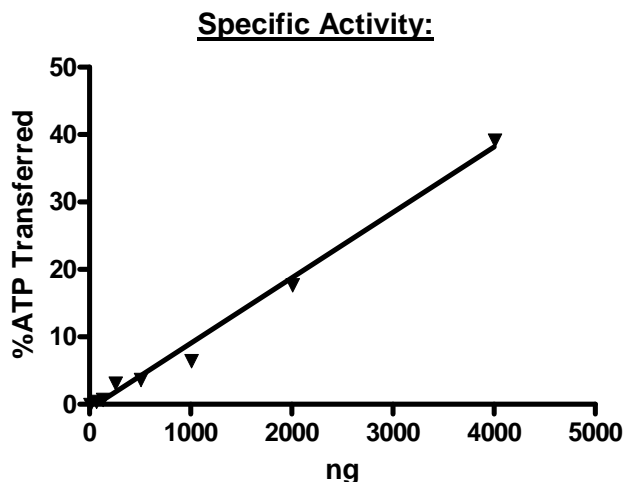
1. Zemann, B. *et al.*, *Blood*. **107(4)**:1454-8 (2006)
2. Maceyka, M. *et al.*, *J Biol Chem*. **280(44)**:37118-29 (2005)

Description: Mouse SPHK2, also known as Sphingosine kinase 2 and SPK2, GenBank Accession No. NM_203280, a.a. 1-617 (full length) with N-terminal His tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=70 kDa.

Specific Activity: 2 pmol/min/ μg
Assay Conditions: Assay was done in Kinase buffer containing 1 mM DTT using 20 μM Sphingosine and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

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

Quality Assurance

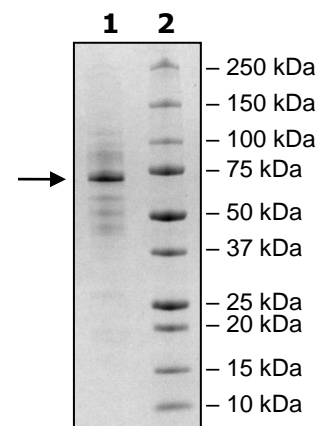


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg SphK2
Lane 2:
Protein Marker

MW: 70 kDa

Purity: $\geq 60\%$



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

Please visit our website at: www.bpsbioscience.com