



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

SPHK1

human, recombinant, , N-terminal HIS tag
Catalog #: 40610

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

Stability: >6 months at -80°C

References:

1. Delon, C., et al., J. Biol. Chem 279 (43) 44763-44774 (2004)
2. Taha, TA et al., FASEB J. 20(3):482-4 (2005)

Description: Human Sphingosine kinase 1 (GenBank Accession No. AF200328), full length with N-terminal His tag, MW=46.9 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. Assay buffer: 50 mM HEPES, pH 7.4, 150 mM NaCl, 5 mM MgCl₂, 1 mM DTT, 3 μM Na-orthovanadate, 0.5 mM ATP, 4 μM D-erythro-sphingosine, and 0.1-0.5 μg/ml sphingosine kinase 1.

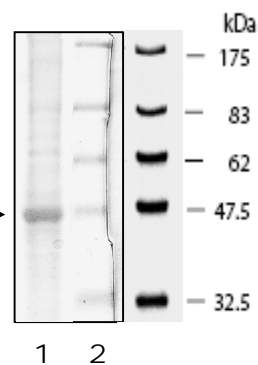
Application:

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

Quality Assurance

10% SDS-PAGE
Coomassie staining

Lane 1:
10 μg SPHK1 →
Lane 2:
Protein Marker
BioLabs (#P7708L)



MW: 46.9 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 **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com