



Echelon Biosciences Inc.
675 Arapeen Drive, Ste 302
Salt Lake City, UT 84108
Phone: 801-588-0455
Fax: 801-588-0497
Email: echelon@echelon-inc.com
<http://www.echelon-inc.com>

Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Sphingosine Kinase 2, active

(Recombinant full-length human protein expressed in Sf9 cells)

Catalog # **E-K069**

Size: 20 µg

Product Description

Sphingosine kinase 2 (sphingosine kinase type 2 isoform, SphK2) is a lipid kinase. Recombinant full-length human sphingosine kinase 2 with N-terminal 6xHis tag was expressed in a baculovirus infected Sf9 insect cell expression system. The GenBank Accession number is AF245447. MW = 69.5 kDa.

Please refer to Certificate of Analysis (COA) for lot specific information about protein concentration, formulation, specific activity, and purity for a certain lot of sphingosine kinase 2 enzyme.

Suggested Assay Condition

Use in Sphingosine Kinase Inhibitor Screen Assay (cat# K-4400): 5 µM D-erythro-sphingosine substrate (cat# K-4401), 5 µM ATP, and 0.05 Unit/mL of recombinant sphingosine kinase 2 in reaction buffer (cat# K-4402 as 10x reaction buffer) for 30 min at room temperature or 37 °C.

Storage and Stability

Store product frozen at or below -70 °C. Enzyme will be stable for at least 6 months at -70 °C as undiluted stock. Flash freeze in working aliquots to avoid multiple freeze/thaw cycles.

Application

Sphingosine kinase 2 is useful for the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

References

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