

Echelon Biosciences Inc.

Sphingosine 1-Phosphate Lyase, active

E-5000P

Support: echelon@echelon-inc.com

Description:

Truncated human sphingosine-1-phosphate lyase (a.a. 62-5686) with N-terminal 6XHis-tag expressed in Sf9 cells and purified using nickel-NTA chromatography.

Applications:

Can be used as a positive control for SPL activity assay, SPL inhibitor screen, & western blot.

Properties:

Size – 2.5 µg & 25 µg

Storage – Store enzyme at -80 °C. Avoid repeated freeze-thaw cycle

Formulation – 10 mM Tris-HCl, 250 mM NaCl, 0.5 mM TCEP, 2.5% glycerol, 1.5 mM methionine, 0.05 mM pyridoxal phosphate, pH 7.4 with proprietary protein stabilizer

Concentration – see certificate of analysis (COA) for lot specific information

Purity – see COA for lot specific information

Activity – see COA for lot specific information

References:

1. Weiler, S.; Braendlin, N.; Beerli, C.; Bergsdorf, C.; Schubart, A.; Srinivas, H.; Oberhauser, B.; Billich, A., Orally active 7-substituted (4-benzylphthalazin-1-yl)-2-methylpiperazin-1-yl]nicotinonitriles as active-site inhibitors of sphingosine 1-phosphate lyase for the treatment of multiple sclerosis. J Med Chem 2014, 57 (12), 5074-84.

Related Products:

| Products | Catalog Number |
|---------------------------|--------------------------------|
| Reagents | |
| Sphingosine 1-Phosphate | S-2000, S-200B, S-200F, S-200T |
| SPL Fluorogenic Substrate | S-200U |
| S1P Beads | S-6110, S-6100 |
| SK I Inhibitor | B-0026 |
| Antibody & Enzyme | |
| S1P Antibody | Z-P300 |
| SPL Overexpressed Lysate | E-5000L |
| Sphingosine Kinase | E-K069, E-K068 |
| Assays | |
| SK Activity Assay | K-3500 |
| S1P ELISA Kit | K-1900 |

Background:

Sphingosine 1-phosphate lyase – SPL (E.C. 4.1.2.27) works as a key regulator in the final step of the sphingolipid degradative pathway by irreversibly breaking down the sphingosine-1-phosphate (S1P) to ethanolamine phosphate and hexadecenal. S1P signaling, through its five receptors, is involved in many biological events such as proliferation, inflammation and apoptosis in cancer and other degenerative diseases.

