
SPHK1

Product code: 11-105

Full-length human SPHK1 [1-384(end) amino acids of accession number NP_001136074.1] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system.

Assay platform : ADP-Glo_Lipid

Substrate : Sphingosine

ATP (μ M) Km app / Bin : 5.9 / 5

Metal : Mg

Reference compound : PF-543

IC50 at ATP Bin (nM) : 17

IC50 at 1 mM ATP (nM) : n.a.
