
SPHK2

Product code: 11-106

Full-length human SPHK2 [1-618(end) amino acids of accession number NP_001191089.1] was expressed as N-terminal GST-fusion protein (92 kDa) using baculovirus expression system.

Assay platform : ADP-Glo_Lipid

Substrate : Sphingosine

ATP (μ M) Km app / Bin : 200 / 200

Metal : Mg

Reference compound : PF-543

IC50 at ATP Bin (nM) : 1400

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