SPHK1

Product code: 11-105

Full-length human SPHK1 [1-384(end) amino acids of accession

fusion protein (69 kDa) using baculovirus expression system.

number NP 001136074.1] was expressed as N-terminal GST-

Reference compound

IC50 at ATP Bin (nM)

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

Assay platform

: ADP-Glo

: Mg

: 17

: PF-543

: Sphingosine

ATP (μ M) Km app / Bin : 5.9 / 5

Metal

Substrate