

---

# ***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. GST-SPHK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Sphingosine

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 20 / 25

Metal : Mg

Reference compound : PF-543

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 3.9

IC<sub>50</sub> at 1 mM ATP (nM) : n.a.

---