

---

# 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
Substrate	: Sphingosine
ATP ( $\mu$ 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.