

---

# EPHA8

Product code: 08-127

Human EPHA8, catalytic domain [571-924 amino acids of accession number NP\_065387.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP ( $\mu$ M) Km app / Bin	: 69 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 61
IC50 at 1 mM ATP (nM)	: 240

