

---

# EPHB4

Product code: 08-131

Human EPHB4, cytoplasmic domain [577-987(end) amino acids of accession number NP\_004435.3] was expressed as N-terminal GST-protein (73 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP ( $\mu$ M) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 230
IC50 at 1 mM ATP (nM)	: 1500

