

---

---

# MER(MERTK)

Product code: 08-108

Human MER, cytoplasmic domain [528-999(end) amino acids of accession number NP\_006334.2] was expressed as N-terminal GST-fusion protein (80 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 36 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.61
IC50 at 1 mM ATP (nM)	: 5.3

