

---

# 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. GST-MER was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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

---