

---

# TRKA(NTRK1)

Product code: 08-186

Human TRKA, cytoplasmic domain [436-790(end) amino acids of accession number NP\_001012331.1] was expressed as N- terminal GST-fusion protein (67 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 65 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.34
IC50 at 1 mM ATP (nM)	: 0.64

---