

---

---

# 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

---

---