

---

---

## TRKB(NTRK2)

Product code: 08-187

Human TRKB, cytoplasmic domain [456-822(end) amino acids of accession number NP\_001018074.1] was expressed as N- terminal GST-fusion protein (69 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 80 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.29
IC50 at 1 mM ATP (nM)	: 0.55

