

---

## TRKC(NTRK3)

Product code: 08-197

Human TRKC, cytoplasmic domain [456-825(end) amino acids of accession number NP\_002521.2] 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	: 47 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.32
IC50 at 1 mM ATP (nM)	: 1.0

