
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. GST-TRKA was purified by using glutathione sepharose chromatography

Assay platform : Mobility Shift Assay
Substrate : CSKtide

ATP (μ M) K_{mapp} / B_{in} : 65 / 75
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 0.34
IC₅₀ at 1 mM ATP (nM) : 0.64
