
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. GST-TRKB was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Src tide

ATP (μ M) K_{mapp} / B_{in} : 80 / 75
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 0.29
IC₅₀ at 1 mM ATP (nM) : 0.55
