
TBK1

Product code 05-115

Full-length human TBK1 [1-729(end) amino acids of accession number NP_037386.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-TBK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CKtide

ATP (μ M) Km app / Bin : 21 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 1.2

IC50 at 1 mM ATP (nM) : 6.2
