

---

# 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.

Assay platform	: MSA
Substrate	: CKtide
ATP ( $\mu$ 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

