

---

# 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

---