

---

# TAK1-TAB1(MAP3K7)

Product code: 09-019

Fused gene of human TAK1 [1-303 amino acids of accession number NP\_663304.1] and human TAB1 [437-504 amino acids of accession number NP\_006107.1] was expressed as N-terminal His-tagged protein (45kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: LRRKtide
ATP ( $\mu$ M) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: 27

