

---

# IKK $\beta$ (IKBK $\beta$ )

Product code: 05-084

Truncated human IKK $\beta$ [1-662 amino acids of accession number NP\_001547.1] was expressed as N-terminal His-tagged protein (77 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Modified I $\kappa$ B $\alpha$ -derived peptide
ATP ( $\mu$ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 410
IC50 at 1 mM ATP (nM)	: >10000