
IKK β (IKBK β)

Product code: 05-084

Truncated human IKK β [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 κ B α -derived peptide
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 410
IC50 at 1 mM ATP (nM)	: >10000
