
IKK ϵ (IKBKE)

Product code: 05-114

Full-length human IKK ϵ [1-716(end) amino acids of accession number NP_054721.1] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: I κ B α peptide
ATP (μ M) Km app / Bin	: 9.5 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.79
IC50 at 1 mM ATP (nM)	: n.a.
