
IKK α (CHUK)

Product code: 05-112

Full-length human IKK α [1-745(end) amino acids of accession number NP_001269.3] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system.

Assay platform	:	IMAP
Substrate	:	I κ B α peptide
ATP (μ M) Km app / Bin	:	41 / 40
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at ATP Bin (nM)	:	310
IC50 at 1 mM ATP (nM)	:	n.a.

