
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.

