
IRAK4

Product code: 09-145

Full-length human IRAK4 [1-460(end) amino acids of accession number NP_057207.2] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: IRAK1 peptide
ATP (μ M) Km app / Bin	: 920 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 11
IC50 at 1 mM ATP (nM)	: 11
