
IRAK1

Product code: 09-101

Truncated human IRAK1 [194-712(end) amino acids of accession number NP_001560.2] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system.

Assay platform	:	IMAP
Substrate	:	SRPKtide
ATP (μ M) Km app / Bin	:	27 / 25
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at ATP Bin (nM)	:	54
IC50 at 1 mM ATP (nM)	:	n.a.

