
PKR(EIF2AK2)

Product code: 05-156

Human PKR , catalytic domain [252-551(end) amino acids of accession number NP_002750.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system.

Assay platform	: IMAP
Substrate	: SRPKtide
ATP (μ M) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 87
IC50 at 1 mM ATP (nM)	: n.a.
