
PEK(EIF2AK3)

Product code: 05-155

Human PEK, cytoplasmic domain [536-1116(end) amino acids and A704S of accession number NP_004827.4] was expressed as N-terminal GST-fusion protein (94 kDa) using E.coli 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)	: 3600
IC50 at 1 mM ATP (nM)	: n.a.

