
PASK

Product code: 02-128

Human PASK, catalytic domain [949-1323(end) amino acids of accession number NP_055963.2] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 9.7 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 13
IC50 at 1 mM ATP (nM)	: 190

