
PRKX

Product code: 01-130

Full-length human PRKX [1-358(end) amino acids of accession number NP_005035.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.59
IC50 at 1 mM ATP (nM)	: n.a.

