
PKC_i(PRKCI)

Product code: 01-139

Full-length human PKC_i[1-587(end) amino acids of accession number NP_002731.3] was expressed as N-terminal GST-fusion protein (94 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 13
IC50 at 1 mM ATP (nM)	: n.a.

