
DCAMKL2

Product code: 02-140

Full-length human DCAMKL2 [1-695(end) amino acids of accession number NP_689832.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 120 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 22
IC50 at 1 mM ATP (nM)	: n.a.

