
MARK3

Product code: 02-122

Full-length human MARK3 [1-729(end) amino acids of accession number NP_002367.5] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.20
IC50 at 1 mM ATP (nM)	: n.a.

