
MARK1

Product code: 02-120

Full-length human MARK1 [1-795(end) amino acids and K794M of accession number NP_061120.3] was expressed as N-terminal GST-fusion protein (116 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.17
IC50 at 1 mM ATP (nM)	: n.a.

