
MRCK α (CDC42BPA)

Product code: 01-107

Truncated human MRCK α [1-574 amino acids of accession number NP_003598.2] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DAPK1tide
ATP (μ M) Km app / Bin	: 0.45 / 1
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: n.a.
