
skMLCK(MYLK2)

Product code: 02-150

Full-length human skMLCK [1-596(end) amino acids of accession number NP_149109.1] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: MLCtide
ATP (μ M) Km app / Bin	: 820 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 51
IC50 at 1 mM ATP (nM)	: 51
