
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
