
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 |
