
CLK2

Product code: 04-127

Full-length human CLK2 [1-499(end) amino acids of accession number NP_001281267.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : DYRKtide-F

ATP (μ M) Km app / Bin : 140 / 150
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 4.6
IC50 at 1 mM ATP (nM) : 28
