
CLK2

Product code 04-127

Full-length human CLK2 [1-499(end) amino acids of accession number AAH53603.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-CLK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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
