
CLK1

Product code: 04-126

Human CLK1, catalytic domain [129-484(end) amino acids of accession number NP_004062.2] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : DYRKtide-F

ATP (μ M) Km app / Bin : 11 / 10
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 9.6
IC50 at 1 mM ATP (nM) : 60
