

---

# 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 ( $\mu$ 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

---