

---

# 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. GST-CLK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : DYRKtide-F

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 11 / 10  
Metal : Mg  
Reference compound : Staurosporine  
IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 9.6  
IC<sub>50</sub> at 1 mM ATP (nM) : 60

---