

---

# 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 ( $\mu$ 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

