

---

# CLK3

Product code: 04-128

Full-length human CLK3 [1-490(end) amino acids of accession number NP\_003983.2] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP ( $\mu$ M) Km app / Bin	: 75 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 820
IC50 at 1 mM ATP (nM)	: n.a.

