

---

# NuaK2

Product code: 02-127

Full-length human NuaK2 [1-628(end) amino acids of accession number NP\_112214.3] was expressed as N-terminal GST-fusion protein (98kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP ( $\mu$ M) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 2.3

