

---

# NuaK1

Product code: 02-126

Full-length human NuaK1 [1-661(end) amino acids of accession number NP\_055655.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CHKtide  
  
ATP ( $\mu$ M) Km app / Bin : 59 / 50  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 1.0  
IC50 at 1 mM ATP (nM) : 3.4

---