

---

# MARK3

Product code 02-122

Full-length human MARK3 [1-729(end) amino acids of accession number NP\_002367.4] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system. GST-MARK3 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform : Mobility Shift Assay  
Substrate : CHKtide

ATP ( $\mu$ M) Km app / Bin : 5 / 5  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.20  
IC50 at 1 mM ATP (nM) : n.a.

---