

---

# MARK4

Product code: 02-123

Full-length human MARK4 [1-688(end) amino acids of accession number NP\_113605.2] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CHKtide  
  
ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 12 / 10  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP  $B_{in}$  (nM) : 0.12  
IC50 at 1 mM ATP (nM) : 5.6

---