

---

# MARK2

Product code: 02-121

Full-length human MARK2 [1-745(end) amino acids of accession number NP\_059672.2] was expressed as N-terminal GST-fusion protein (110 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP ( $\mu$ M) Km app / Bin	: 8.8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.12
IC50 at 1 mM ATP (nM)	: n.a.

