

---

---

# MELK

Product code: 02-124

Truncated human MELK [1-493 amino acids of accession number NP\_055606.1] was expressed as N-terminal GST-fusion protein (83 kDa) using E. coli expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP ( $\mu$ M) Km app / Bin	: 38 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.81
IC50 at 1 mM ATP (nM)	: n.a.

