

---

---

# MRCK $\beta$ (CDC42BPB)

Product code: 01-108

Truncated human MRCK $\beta$  [1-473 amino acids of accession number NP\_006026.3] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DAPK1tide
ATP ( $\mu$ M) Km app / Bin	: 0.67 / 1
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.6
IC50 at 1 mM ATP (nM)	: n.a.

---

---