

---

# 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.

---