

---

# ROCK2

Product code: 01-110

Human ROCK2, catalytic domain [1-553 amino acids of accession number NP\_004841.2] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : LIMKtide  
  
ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 7.4 / 5  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP  $B_{in}$  (nM) : 0.92  
IC50 at 1 mM ATP (nM) : 21

---