

---

# DAPK1

Product code: 02-134

Human DAPK1, catalytic domain [1-289 amino acids of accession number NP\_004929.2] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : DAPK1tide  
  
ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 1.1 / 1  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP  $B_{in}$  (nM) : 2.9  
IC50 at 1 mM ATP (nM) : 490

---