

---

---

# HPK1(MAP4K1)

Product code: 07-410

Human HPK1, catalytic domain [1-346 amino acids of accession number NP\_009112.1] was expressed as N-terminal DYKDDDDK tagged protein (41 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : S6K2 peptide  
  
ATP ( $\mu$ M) Km app / Bin : 22 / 25  
Metal : Mg  
Reference compound : K252a  
IC50 at ATP Bin (nM) : 6.9  
IC50 at 1 mM ATP (nM) : 32

---

---