

---

# WNK3

Product code: 05-181

Human WNK3, catalytic domain [1-434 amino acids of accession number NP\_065973.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : SPAKtide  
  
ATP ( $\mu$ M)  $K_m$  app / Bin : 48 / 50  
Metal : Mg+Mn  
Reference compound : K252a  
IC50 at ATP Bin (nM) : 1300  
IC50 at 1 mM ATP (nM) : n.a.

---