

---

# WNK2

Product code: 05-180

Human WNK2, catalytic domain [166-489 amino acids of accession number NP\_006639.3] was expressed as N-terminal GST-fusion protein (65 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) : 2300  
IC50 at 1 mM ATP (nM) : n.a.

---