

---

# WNK1

Product code: 05-179

Human WNK1, catalytic domain [1-491 amino acids of accession number NP\_061852.3] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: SPAKtide
ATP ( $\mu$ M) $K_m$ app / Bin	: 140 / 150
Metal	: Mg+Mn
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 4500
IC50 at 1 mM ATP (nM)	: n.a.

