

---

# 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) Km app / Bin	: 48 / 50
Metal	: Mg+Mn
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 1300
IC50 at 1 mM ATP (nM)	: n.a.

