

---

# 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) Km app / Bin	: 48 / 50
Metal	: Mg+Mn
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 2300
IC50 at 1 mM ATP (nM)	: n.a.

