

---

# TNIK

Product code: 07-138

Human TNIK, catalytic domain [1-314 amino acids of accession number NP\_055843.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Moesin-derived peptide
ATP ( $\mu$ M) $K_m$ app / $B_{in}$	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP $B_{in}$ (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 11

---