
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 (μ M) K_m app / B_{in}	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP B_{in} (nM)	: 1.0
IC ₅₀ at 1 mM ATP (nM)	: 11
