
TXK

Product code: 08-183

Human TXK, catalytic domain [260-527(end) amino acids of accession number NP_003319.2] was expressed as N-terminal GST-fusion protein (58 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 110 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 45
IC50 at 1 mM ATP (nM)	: 220

