
TEC

Product code: 08-182

Human TEC, catalytic domain [359-631(end) amino acids of accession number NP_003206.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 55 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 20
IC50 at 1 mM ATP (nM)	: 220

