
ITK

Product code: 08-181

Full-length human ITK [2-620(end) amino acids of accession number NP_005537.3] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 6.1 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.2
IC50 at 1 mM ATP (nM)	: 200

