
BTK

Product code: 08-180

Full-length human BTK [2-659(end) amino acids of accession number NP_000052.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 22 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 24
IC50 at 1 mM ATP (nM)	: 93

