
ZAP70

Product code: 08-177

Full-length human ZAP70 [1-619(end) amino acids of accession number NP_001070.2] was expressed as N-terminal GST-fusion protein (97 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 3.3 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.76
IC50 at 1 mM ATP (nM)	: 34
