
FYN[isoform b]

Product code: 08-531

Full-length human FYN [isoform b] [1-534(end) amino acids of accession number NP_694592.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.8
IC50 at 1 mM ATP (nM)	: 42

