
FYN[isoform a]

Product code: 08-168

Full-length human FYN [isoform a] [1-537(end) amino acids of accession number NP_002028.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 36 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.9
IC50 at 1 mM ATP (nM)	: 24

