
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 : Src tide

ATP (μ M) K_m app / B_{in} : 20 / 25

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 2.8

IC₅₀ at 1 mM ATP (nM) : 42
