
SYK

Product code 08-176

Full-length human SYK [1-635(end) amino acids of accession number NP_003168.2] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system. GST-SYK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 59 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.29

IC50 at 1 mM ATP (nM) : 0.63
