
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.

Assay platform : MSA
Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 59 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.4
IC50 at 1 mM ATP (nM) : 4.5
