

---

# 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 ( $\mu$ 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

