

---

# LYNb

Product code 08-172

Full-length human LYNb [1-491(end) amino acids of accession number NP\_001104567.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-LYNb was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform : Mobility Shift Assay

Substrate : Srctide

ATP ( $\mu$ M) Km app / Bin : 18 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.0

IC50 at 1 mM ATP (nM) : 21

---