

---

# LYNa

Product code: 08-171

Full-length human LYNa [1-512(end) amino acids of accession number NP\_002341.1] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Srctide

ATP ( $\mu$ M) Km app / Bin : 14 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.0

IC50 at 1 mM ATP (nM) : 22

---