
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.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : Srctide |
| ATP (μ 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 |

