
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. GST-LYNa was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Srctide

ATP (μ M) Kmapp / Bin : 14 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.0

IC50 at 1 mM ATP (nM) : 22
