
PIM2

Product code 02-155

Full-length human PIM2 [1-311(end) amino acids of accession number NP_006866.2] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system. GST-PIM2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : S6K2 peptide

ATP (μ M) Km app / Bin : 4 / 5

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 14

IC50 at 1 mM ATP (nM) : 480
