
PIM1

Product code: 02-054

Full-length human PIM1 [1-313(end) amino acids of accession number NP_002639.1] was expressed as N-terminal His-tagged protein (39 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : S6K2 peptide

ATP (μ M) K_m app / B_{in} : 640 / 500
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP B_{in} (nM) : 10
IC50 at 1 mM ATP (nM) : 20
