
PIM3

Product code: 02-156

Full-length human PIM3 [1-326(end) amino acids and V300A of accession number NP_001001852.2] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : S6K2 peptide

ATP (μ M) K_m app / B_{in} : 130 / 150
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 0.36
IC₅₀ at 1 mM ATP (nM) : 0.71
