

---

---

# 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 ( $\mu$ M) Km app / Bin	: 640 / 500
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 20

