

---

# 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.

Assay platform	: MSA
Substrate	: S6K2 peptide
ATP ( $\mu$ M) Km app / Bin	: 4 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 14
IC50 at 1 mM ATP (nM)	: 480

---