

---

---

## PYK2(PTK2B)

Product code: 08-138

Full-length human PYK2 [1-967(end) amino acids of accession number NP\_775267.1] was expressed as N-terminal GST-fusion protein (138 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP ( $\mu$ M) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 4.9

