

---

# TYK2

Product code: 08-147

Human TYK2, catalytic domain [871-1187(end) amino acids of accession number NP\_003322.3] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 18 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 7.0

---