

---

# JAK3

Product code: 08-046

Human JAK3, catalytic domain [795-1124(end) amino acids of accession number NP\_000206.2] was expressed as N-terminal His-tagged protein (41 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 3.5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.20
IC50 at 1 mM ATP (nM)	: 12

