

---

# JAK2

Product code: 08-045

Human JAK2, catalytic domain [826-1132(end) amino acids of accession number NP\_004963.1] was expressed as N-terminal His-tagged protein (39 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : Srctide  
  
ATP ( $\mu$ M) Km app / Bin : 13 / 10  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.34  
IC50 at 1 mM ATP (nM) : 6.0

---