

---

# JAK1

Product code: 08-144

Human JAK1, catalytic domain [850-1154(end) amino acids of accession number NP\_002218.2] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : JAK1 substrate peptide  
  
ATP ( $\mu$ M) Km app / Bin : 68 / 75  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.71  
IC50 at 1 mM ATP (nM) : 10

---