

---

# 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. His-tagged JAK3 was purified by using Ni-NTA affinity chromatography and gel filtration chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP ( $\mu\text{M}$ )  $K_{\text{mapp}}$  /  $B_{\text{in}}$  : 3.5 / 5

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{\text{in}}$  (nM) : 0.20

IC<sub>50</sub> at 1 mM ATP (nM) : 12

---