
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 (μ 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
