
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 : Src tide

ATP (μM) K_m app / B_{in} : 13 / 10

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.34

IC₅₀ at 1 mM ATP (nM) : 6.0
