
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.

Assay platform : MSA
Substrate : Srctide
ATP (μ M) Km app / Bin : 3.5 / 5
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 0.20
IC50 at 1 mM ATP (nM) : 12
