
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 (μ 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
