

PRODUCT DATA SHEET



JAK2

CATALOG NO.: SE-435

LOT NO.: Template

DESCRIPTION: MW=42.5 kDa; human recombinant JAK2 (GenBank Accession No. NM_004972), (a.a. 808-end) with N-terminal His tag, expressed in a Baculovirus infected Sf9 cell expression system

PURITY: >80% by SDS-PAGE.

SPECIFIC ACTIVITY: 4000 U/mg. One unit is defined as the amount of enzyme that will phosphorylate 1 nmol of tyrosine substrate per minute at pH 7.4 and 30°C. Assay buffer: 50 mM HEPES, pH 7.4, 3 mM MgCl₂, 3 mM MnCl₂, 1 mM DTT, 3 μM Na-orthovanadate, 0.5 mM ATP, 30 μg/ml Poly (Glu:Tyr)4:1 substrate, and 4 μg/ml recombinant JAK2.

APPLICATIONS: Study enzyme regulation and kinetics, phosphorylate target substrates, screen for inhibitors.

SUPPLIED AS: 10 μg at XXX mg/ml in 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% glycerol, 3 mM DTT.

STORAGE: -80°C. After initial defrost, aliquot product into individual tubes and refreeze the remaining, unused enzyme quickly by snap-freezing in a dry/ice ethanol bath or liquid nitrogen, if possible. Avoid repeated freeze/defrost cycles.

NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.