

STAT1 beta Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # S52-54G

Lot # S203-1

Product Description

Recombinant full-length human STAT1β was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_139266](#).

Gene Aliases

ISGF-3; STAT91; DKFZp686B04100

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

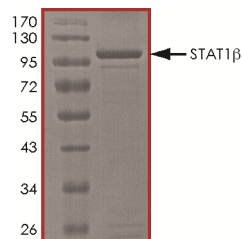
Scientific Background

STAT1β is a member of the signal transducers and activators of transcription (STAT) family of proteins that carry out a dual function: signal transduction and activation of transcription. STAT1β transcription factor is specific for the IFN pathway and plays a central role in mediating many, if not all, IFN-dependent biological responses (1). Presence of STAT1β leads to an efficient antiviral response when cells were infected with virus suggesting that a STAT-dependent pathway is activated following virus infection by endogenously produced IFN. Virus-induced STAT protein translocation from the cytoplasmic compartment can be detected within 3 h of infection.

References

1. Improti, T. et al: Susceptibility to virus infection is determined by a Stat-mediated response to the autocrine effect of virus-induced type I interferon. Cytokine. 1997 Jun;9(6):383-93.

Purity



The purity of STAT1 beta was determined to be **>90%** by densitometry. Approx. MW **118kDa**.

Upstream Active Kinases

SRC, Active
MEK1, Active

Cat # S19-10G
Cat # M02-13G

STAT1 beta Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	S52-54G
Specific Lot Number	S203-1
Purity	>90%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.