

Catalogue #	Aliquot Size
<b>S52-54G-20</b>	<b>20 µg</b>
<b>S52-54G-50</b>	<b>50 µg</b>

## STAT1 beta Protein

Full-length recombinant protein expressed in Sf9 cells

### Catalog # S52-54G

Lot # S131-1

### Product Description

Recombinant full-length human STAT1 $\beta$  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.

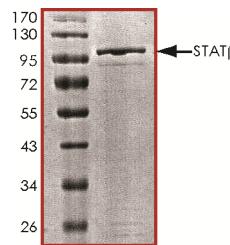
### Scientific Background

STAT1 $\beta$  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. STAT1 $\beta$  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 STAT1 $\beta$  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. Impronta, 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 was determined to be **>85%** 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	S131-1
Purity	>85%
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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