

STAT3

Full length recombinant protein expressed in Sf9 cells

Catalog # S54-54G

Lot # G323-1

Product Description

Recombinant full-length human STAT3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_139276](#).

Gene Aliases

APRF, FLJ20882, MGC16063

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 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.

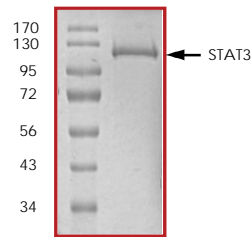
Scientific Background

STAT3 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. STAT3 is widely expressed and becomes activated through phosphorylation on tyrosine as a DNA binding protein in response to a variety of stimuli such as EGF, IL-6, PDGF, IL-2 and G-CSF (1). This phosphoprotein forms homodimers as well as heterodimers with other members of the STAT family and translocate to the nucleus in order to modulate the transcription of various genes (2).

References

1. Zhong, Z. et al: Stat3: a STAT family member activated by tyrosine phosphorylation in response to epidermal growth factor and interleukin-6. *Science*. 1994 Apr 1;264(5155):95-8.
2. Tian, S. S. et al: Rapid activation of the STAT3 transcription factor by granulocyte colony-stimulating factor. *Blood*. 1994 Sep 15;84(6):1760

Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **120kDa**.

Upstream Active Kinases

SRC, Active
MEK1, Active

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

STAT3

Full length recombinant protein expressed in Sf9 cells

Catalog Number

S54-54G

Specific Lot Number

G323-1

Purity

>90%

Concentration

0.2 $\mu\text{g}/\mu\text{l}$

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.