

STAT5 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # S56-54H

Lot # S119-1

Product Description

Recombinant full-length human STAT5 was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_003152](#).

Gene Aliases

MGF, STAT5A

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 2mM DTT, 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

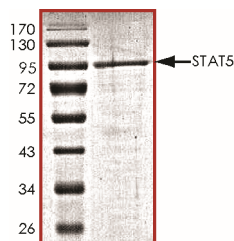
STAT5 (signal transducers and activators of transcription 5) is a member of the STAT family and this protein is phosphorylated by receptor associated protein kinases in response to cytokines and growth factors (1). The differentiation of T helper (Th) cells is regulated by members of the STAT family of signaling molecules. In BCR-ABL-positive cells, STAT5 is constitutively activated by tyrosine phosphorylation. STAT5 activation results in upregulation of Bcl-X_L and increased resistance to induction of apoptosis. STAT5 is involved in expression and growth hormone-mediated sexually dimorphic regulation of cytochrome P450 3A10/lithocholic acid 6 beta-hydroxylase (2).

References

1. Wang, D. et al: Naturally occurring dominant negative variants of Stat5. *Molec. Cell Biol.* 16: 6141-6148, 1996.
2. A Subramanian et al: STAT 5 and NF-Y are involved in expression and growth hormone-mediated sexually dimorphic regulation of cytochrome P450 3A10/lithocholic acid 6beta-hydroxylase. *Nucleic Acids Res.* 1998; 26(9): 2173-2178.

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

Purity



The purity of STAT5 was determined to be **>90%** by densitometry. Approx. MW **92kDa**.

STAT5 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	S56-54H
Lot #	S119-1
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.