

STAT4 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # S55-54H

Lot # Q175-3

Product Description

Recombinant full-length human STAT4 was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_003151](#).

Gene Aliases

None

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM 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

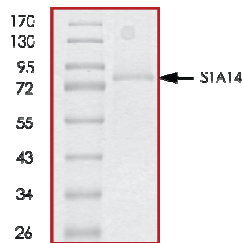
STAT4 (signal transducer and activator of transcription 4) is a member of the STAT family that is phosphorylated by receptor-associated protein kinases. STATs form homo- or heterodimers and translocate to the cell nucleus where they act as transcription activators. T helper (Th) cells differentiation is regulated by STAT4 (1). STAT4 controls T1/T2 balance during the response to live influenza virus and isolated viral proteins. The differentiation of gamma interferon (IFN-γ)-producing Th1 and Tc1 cells after inoculation of live virus occurred independently of STAT4 expression. STAT4 has been shown to control pathogenesis of chronic obstructive pulmonary disease (2).

References

1. Andrea L W. et al: The biology of Stat4 and Stat6. *Oncogene*, 2000,19: 21, 2577-2584.
2. Cosio. G M.: Autoimmunity, T-cells and STAT-4 in the pathogenesis of chronic obstructive pulmonary disease. *Eur Respir J*, 2004; 24:3-5.

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 STAT4 was determined to be **>95%** by densitometry. Approx. MW **87kDa**.

STAT4 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	S55-54H
Specific Lot Number	Q175-3
Purity	>95%
Concentration	0.2 µg/µl
Stability	1 yr 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.