

PRODUCT INFORMATION



STAT1 β (human, recombinant)

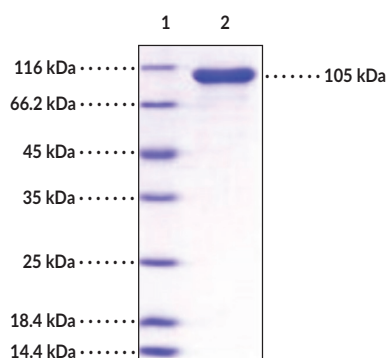
Item No. 32022

Overview and Properties

Synonyms:	Signal Transducer and Activator of Transcription-1 β , STAT84, Transcription Factor ISGF-3 Component p84
Source:	Recombinant human N-terminal His- and GST-tagged STAT1 β expressed in insect cells (baculovirus)
Amino Acids:	1-712 (full length)
Molecular Weight:	111 kDa
Storage:	-80°C (as supplied)
Stability:	≥ 1 year
Purity:	$\geq 92\%$ estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 20 mM Tris, pH 7.4, with 150 mM sodium chloride
Bioactivity:	See figures for details

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: STAT1 β

SDS-PAGE analysis of STAT1 β . This protein has a calculated molecular weight of 111 kDa. It has an apparent molecular weight of approximately 105 kDa by SDS-PAGE under reducing conditions.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/15/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

STAT1, also known as STAT1 α , is a transcription factor and member of the STAT protein family with roles in innate and adaptive immunity.¹ It is composed of an N-terminal domain that is essential to protein-protein interactions and dimerization, a DNA binding domain that facilitates nuclear import and export, as well as DNA binding, a linker domain, tail segment, and a transactivation domain that facilitates transcription of target genes.² STAT1 β is an isoform of STAT1 that is formed by alternative splicing and lacks the 38-amino acid transactivation domain.³ Upon phosphorylation by JAKs, STAT1 dimerizes and is translocated to the nucleus to activate transcription of IFN- γ -inducible genes.⁴ Because the C-terminal transactivation domain is required for STAT1 interaction with the transcriptional coactivator CREB-binding protein (CBP)/p300 and gene transcription, STAT1 β is considered a dominant-negative regulator of STAT1 and overexpression of STAT1 β inhibits IFN- γ -induced gene expression in RAW 264.7 cells. STAT1 β phosphorylation is increased and IFN- γ -induced JAK1, JAK2, and STAT1 activation is decreased in RAW 264.6 cells infected with *L. major* and *L. mexicana* parasites, as well as *M. avium* bacteria.^{4,5} STAT1 β protein levels are reduced in patient-derived esophageal squamous cell carcinoma (ESCC) tumor samples and are positively correlated with lymph node metastasis, invasion, and shorter overall survival.³ Cayman's STAT1 β (human, recombinant) protein consists of 950 amino acids and has a calculated molecular weight of 111 kDa.

References

1. Meissl, K., Macho-Maschler, S., and Müller, M. The good and the bad faces of STAT1 in solid tumours. *Cytokine* **89**, 12-20 (2017).
2. Boisson-Dupuis, S., Kong, X.-F., Okada, S., *et al.* Inborn errors of human STAT1: Allelic heterogeneity governs the diversity of immunological and infectious phenotypes. *Curr. Opin. Immunol.* **24(4)**, 364-378 (2012).
3. Zhang, Y., Chen, Y., Yun, H., *et al.* STAT1 β enhances STAT1 function by protecting STAT1 α from degradation in esophageal squamous cell carcinoma. *Cell Death Dis.* **8(10)**, e3077 (2017).
4. Alvarez, G.R., Zwilling, B.S., and Lafuse, W.P. Mycobacterium avium inhibition of IFN- γ signaling in mouse macrophages: Toll-like receptor 2 stimulation increases expression of dominant-negative STAT1 β by mRNA stabilization. *J. Immunol.* **171(12)**, 6766-6773 (2003).
5. Bhardwaj, N., Rosas, L.E., Lafuse, W.P., *et al.* *Leishmania* inhibits STAT1-mediated IFN- γ signaling in macrophages: Increased tyrosine phosphorylation of dominant negative STAT1 β by *Leishmania mexicana*. *Int. J. Parasitol.* **35(1)**, 75-82 (2005).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM