

NFATC1 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # N12-30G

Lot # X609-1

Product Description

Recombinant full-length human NFATC1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_172390](#).

Gene Aliases

NF-ATC, NFATc, NFAT2

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.

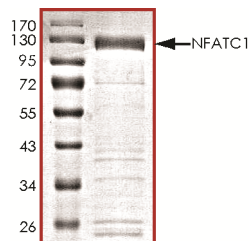
Scientific Background

NFATC1 is a member of the NFAT family of proteins which are Ca²⁺/calcineurin-responsive transcription factors primarily recognized for their central roles in T lymphocyte activation and cardiac valve development (1). NFAT consists of at least two components: a preexisting cytosolic component that translocates to the nucleus upon T-cell receptor (TCR) stimulation, and an inducible nuclear component. Proteins belonging to the NFAT family of transcription factors play a central role in inducible gene transcription during immune response. The product of NFATC1 is an inducible nuclear component and functions as a major molecular target for the immunosuppressive drugs such as cyclosporin A (2).

References

- Buchholz, M. et al: Overexpression of c-myc in pancreatic cancer caused by ectopic activation of NFATc1 and the Ca²⁺/calcineurin signaling pathway. EMBO. J. 2006; 25(15):3714-24.
- Jhun, B.S. et al: Inhibition of AMP-activated protein kinase suppresses IL-2 expression through down-regulation of NF-AT and AP-1 activation in Jurkat T cells. Biochem. Biophys. Res. Commun. 2006; 351(4):986-92.

Purity



The purity of NFATC1 was determined to be **>90%** by densitometry. Approx. MW **125kDa**.

NFATC1 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	N12-30G
Specific Lot Number	X609-1
Purity	>90%
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.

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.