

FOS Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # F66-30G

Lot # Q355-3

Product Description

Recombinant full-length human FOS was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_005252](#).

Gene Aliases

c-FOS

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

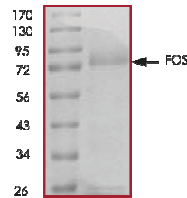
FOS is a member of the FOS gene family that consists of 4 members: FOS, FOSB, FOSL1, and FOSL2. FOS gene encodes a leucine zipper protein that can dimerize with proteins of the JUN family, thereby forming the transcription factor complex AP-1. FOS proteins have been implicated as regulators of cell proliferation, differentiation, and transformation (1). In some cases, expression of the FOS gene has also been associated with apoptotic cell death. FOS overexpression leads to decreased phosphorylation and dimerization of STAT1, which in turn downregulates p21 gene expression (2). This regulatory pathway may enhance the proliferation of lymphocytes in rheumatoid arthritis patients.

References

1. Saez, E. et al: c-fos is required for malignant progression of skin tumors. *Cell* 82: 721-732, 1995.
2. Hikasa, M. et al: p21(waf1/cip1) is down-regulated in conjunction with up-regulation of c-Fos in the lymphocytes of rheumatoid arthritis patients. *Biochem. Biophys. Res. Commun.* 304: 143-147, 2003.

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 FOS was determined to be **>90%** by densitometry. Approx. MW **78kDa**.

Upstream Active Kinases

p38 α
p38 β

Catalog # M39-10G
Catalog # M36-10G

FOS Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number **F66-30G**
Specific Lot Number **Q355-3**

Purity
Concentration
Stability
Storage & Shipping

>90%
0.2 µg/µl
1 yr At -70°C from date of shipment
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.