

Recombinant c-Fos protein

Catalog No: 31115

Expressed In: Baculovirus

Quantity: 5 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: 5 µg of Recombinant c-Fos in Dilution Buffer AM1 (20 mM Tris-Cl (pH 7.5), 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA).

Background: **c-Fos** is one of the proteins that form the heteromeric **AP-1 transcription factor complex**. AP-1 proteins play a role in the expression of many genes involved in the regulation of cellular processes such as differentiation, proliferation and apoptosis. The transcription factor AP-1 is composed of a mixture of heterodimeric protein complexes derived from the Fos and Jun families, including **c-Fos, FosB, Fra-1, c-Jun, JunB and JunD**. c-Fos is a nuclear phosphoprotein that forms a tight but non-covalently linked complex with the Jun/AP-1 transcription factor. In the heterodimer, Fos and Jun/AP-1 basic regions each seem to interact with symmetrical DNA half sites. Upon TGF-β activation, c-Fos forms a multimeric SMAD3/SMAD4/Jun/Fos complex at the AP-1/SMAD-binding site to regulate TGF-β-mediated signaling. c-Fos has a critical function in regulating bone morphogenesis.

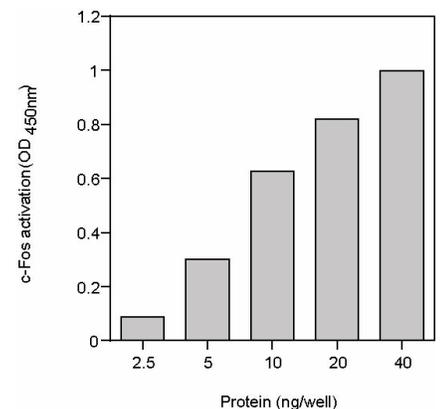
Protein Details: Recombinant c-Fos was expressed from full length (accession number NM 003131) with an N-terminal His-Tag in a baculovirus system and purified by an affinity column in combination with FPLC chromatography. The purified recombinant protein is greater than 90% homogeneous and contains no detectable protease, DNase and RNase activity.

Application Notes: Recombinant c-Fos is suitable for DNA and protein-protein interaction assays. 100 ng is sufficient for DNA-protein and protein-protein interaction studies. The molecular weight of the protein is ~50 kDa. The standard curve for TransAM® AP-1 c-Fos was generated using the range of 40-2.5 ng of protein.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



TransAM® standard curve generated using Recombinant c-Fos protein. The standard curve for TransAM® was generated using a range of 40-10 ng of protein and run on the TransAM® AP-1 c-Fos ELISA Kit.