

Recombinant Human TGF Beta-3 (TGFB3)

Recombinant human protein expressed in E. coli cells

Catalog # T833-40N

Lot # G3223-16

Product Description

Recombinant Human TGF Beta-3 (TGFB3) was expressed in E. coli cells. The protein accession number is [P10600](#)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

The filter sterilized clear protein solution contains 20% Ethanol and 0.12% Acetic acid.

Storage and Stability

This cytokine can be stored in working aliquots at 2° - 8°C for two weeks. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

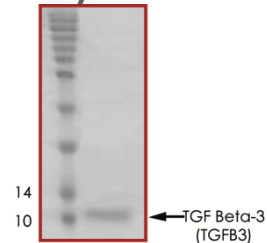
Scientific Background

Transforming Growth Factor Beta-3 is a disulfide-linked homodimeric, non-glycosylated, protein that is one of the isoforms of the highly pleiotropic cytokines that are secreted by all cell types. TGF Beta-1, TGF Beta-2, and TGF Beta-3 use the same receptor and act as cellular switches that regulate cell growth, proliferation, differentiation, motility, extracellular matrix synthesis, and immune function regulation. Defects of this gene can cause arrhythmic right ventricular dysplasia 1 and rienhoff syndrome. TGFB3 first interacts with the accessory receptor betaglycan (TGF Beta-R3) and a type II serine/threonine kinase receptor (TGF Beta-R2). The receptor then activates a type I serine/threonine kinase receptor such as ALK-1 or TGF Beta-R1 (ALK-5) through phosphorylation, and the activated type I receptor activates Smad proteins to regulate transcription.

Specific Activity

The ED50 is 0.01-0.04 ng/mL as determined by its ability to inhibit IL-4 dependent proliferation of HT-2 mouse cells

Purity



Recombinant Human TGF Beta-3 (TGFB3) resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **13.0 kDa**

Recombinant Human TGF Beta-3 (TGFB3)

Recombinant human protein expressed in E. coli cells

Catalog Number **T833-40N**

Specific Lot Number **G3223-16**

Purity **95%**

Stability **6 mos. at -20°C from date of shipment**

Storage & Shipping

This cytokine can be stored in working aliquots at 2° - 8°C for two weeks. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

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.