

TGF- β

Recombinant Human Transforming Growth Factor Beta 1

Catalog # rhtgf-b

For research use only, not for diagnostic or therapeutic use

Version # 10B18-MM

PRODUCT INFORMATION

Content:

- 10 μ g of TGF- β provided as a sterile filtered white lyophilized powder.

Storage and stability:

- TGF- β is shipped at room temperature. Upon receipt it should be stored at -20°C.
- Upon resuspension, TGF- β should be aliquoted and stored at 4°C for short term storage or -20°C for long term storage. Product is stable one week at 4°C and 12 months at -20°C when properly stored. For long term storage, it is recommended to add a carrier protein (0.1% HSA or BSA).

Note: Avoid repeated freeze-thaw cycles.

DESCRIPTION

Tumor growth factor-beta (TGF- β) belongs to a family of structurally related cytokines that regulate a plethora of cellular functions, such as proliferation, apoptosis, differentiation and migration. Recombinant human TGF- β is a 112 amino acid protein. TGF- β is produced in *E. coli*.

Quality control

- Purity greater than 95% as determined by SDS-PAGE

CHARACTERISTICS

Working Concentration: 0.1 - 10 ng/ml

Solubility

It is recommended to reconstitute lyophilized TGF- β in sterile water at a concentration of not less than 100 μ g/ml, which can then be further diluted in aqueous solutions.

TECHNICAL SUPPORT

Toll free (US): 888-457-5873
Outside US: (+1) 858-457-5873
Europe: +33 562-71-69-39
E-mail: info@invivogen.com
Website: www.invivogen.com


3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA