TGF-β3 (mouse), (recombinant)

The transforming growth factors (TGFs) are multifunctional peptides that regulate growth and differentiation in a variety of cells. Recent data suggests that individual TGF- β isoforms (TGF- β 1, - β 2 and - β 3) have overlapping, yet distinct biological actions and target cell specificities, both in developing and adult tissues. TGF- β 3 is a new isoform that is presumed to play an important role in wound repair and scarring. TGF- β 3 is also thought to be involved in osteoblast proliferation, chemotaxis, and collagen synthesis.

- High purity
- Carrier-free

Ordering Information

Order Online »

ALX-201-817-0010	10µg
ALX-201-817-0100	100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable at +4°C for 1 year from date of purchase.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Transforming growth factor β3

Appearance Clear solution.

Biological Activity The activity is determined by the cell toxicity assay, using

the WHO Standard 98/608 as a direct comparison, which

is typically observed at concentations <0.05ng/ml.

Concentration 0.25mg/ml

Endotoxin Content ≤1 EU/μg protein measured by kinetic LAL analysis.

Formulation Liquid. In 20% ethanol and 0.12% acetic acid.

MW ~25.5kDa

Purity ≥98% (Reducing and Non-reducing SDS-PAGE)

Source Produced in E. coli. Non-glycosylated, disulfide-linked

homodimer, containing two 112 amino acid chains.

UniProt ID P17125

Last modified: May 29, 2024

