

TGF- β 3 (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.

Ordering Information

[Order Online »](#)

ALX-201-817-0010	10 μ g
ALX-201-817-0100	100 μ g

Manuals, SDS & CofA

[View Online »](#)

- High purity
- Carrier-free

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 β 3

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 concentrations <0.05ng/ml.

Concentration 0.25mg/ml

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

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

MW ~25.5kDa

Purity $\geq 98\%$ (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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)