

TGF- β 1 (human), (recombinant)

Transforming growth factor β 1 (TGF- β 1) is a member of the TGF- β superfamily of cytokines. Members of this family exhibit regulatory activity in immunity, proliferation, adhesion, migration, and growth inhibition pathways. TGF- β 1 signals through SMAD proteins via the TGF- β RI and TGF- β RII receptors.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT221-0020	20 μ g
-----------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Differentiation inhibiting factor, Cartilage-inducing factor, Transforming growth factor β 1
Endotoxin Content	≤ 1 EU/ μ g protein (LAL test)
Formulation	Lyophilized. Contains 0.1% trifluoroacetic acid and trehalose (in a 20:1 trehalose to protein ratio).
MW	Dimer $\sim 12.8/25.6$ kDa
Purity	$\geq 95\%$ (SDS-PAGE)
Reconstitution	Reconstitute in sterile 10 mM HCl at 0.1 mg/mL.
Source	Produced in CHO cells. Contains 112/224 amino acids.
UniProt ID	P01137



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)