

Growth and differentiation factor-15 (human), (recombinant)

Growth and differentiation factor 15 (GDF-15) is a TGF- β family member, made by the placenta and heart tissues, that has a role in regulating inflammatory and apoptotic pathways. GDF-15 has become an emerging marker of early heart disease and has the potential as being used as a molecule for screening patients for early heart failure.

Ordering Information

[Order Online »](#)

ALX-201-803-0005	5 μ g
------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

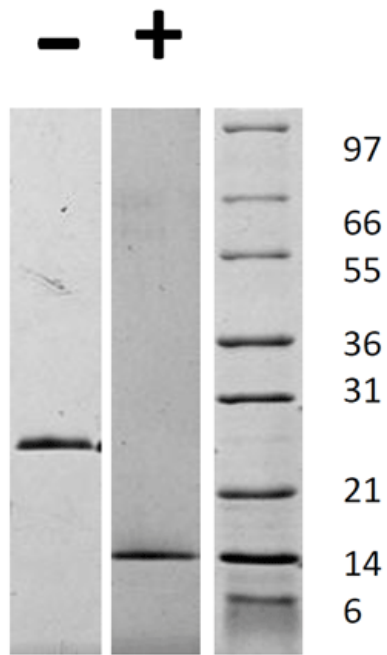


Figure 1: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human GDF-15 is predicted to be a homodimer with a predicted MW of 24.5 kDa.

Handling & Storage

Use/Stability	It is recommended that a carrier protein (0.1% HSA or BSA) is added for long term storage.
Handling	Centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GDF-15, MIC-1, Placental TGF- β , Prostate differentiation factor
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the inhibition of DU-145 cells which is typically observed at a concentration of 1-2 μ g/ml.
Endotoxin Content	≤ 1 EU/ μ g protein measured by kinetic LAL analysis.
Formulation	Lyophilized. Sterile filtered.
MW	~24.5kDa
Purity	$\geq 95\%$ (HPLC, Reducing and Non-reducing SDS-PAGE, UV spectroscopy at 280 nm)
Reconstitution	When reconstituting the product, gently pipet and wash down the sides of the vial to ensure full recovery of the protein into solution. It is recommended to reconstitute the lyophilized product with sterile 5 mM acetic acid (AcOH) at a concentration of 0.1 mg/ml, which can be further diluted into other aqueous solutions.
Source	Produced in <i>E. coli</i> . Non-glycosylated, disulfide linked homodimer, containing two identical 113 amino acid chains.
UniProt ID	Q99988



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)