

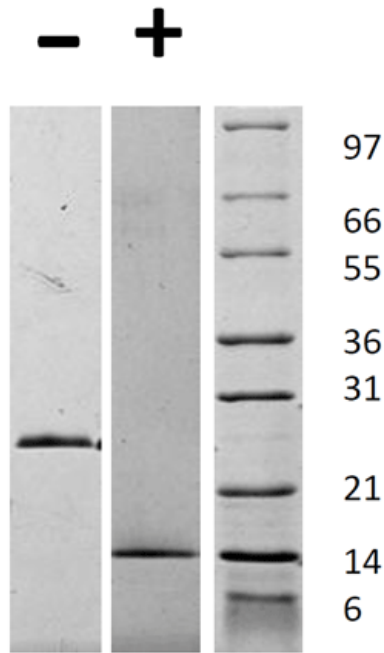
# Growth and differentiation factor-15 (human), (recombinant)

Growth and differentiation factor 15 (GDF-15) is a TGF- $\beta$  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.

**Manuals, SDS & CofA**

[View Online »](#)

- Carrier-free



**Figure 1:** 1 $\mu$ 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

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	GDF-15, MIC-1, Placental TGF- $\beta$ , Prostate differentiation factor
<b>Appearance</b>	White lyophilized (freeze-dried) powder.
<b>Biological Activity</b>	The activity is determined by the inhibition of DU-145 cells which is typically observed at a concentration of 1-2 $\mu$ g/ml.
<b>Endotoxin Content</b>	$\leq$ 1 EU/ $\mu$ g protein measured by kinetic LAL analysis.
<b>Formulation</b>	Lyophilized. Sterile filtered.
<b>MW</b>	~24.5kDa
<b>Purity</b>	$\geq$ 95% (HPLC, Reducing and Non-reducing SDS-PAGE, UV spectroscopy at 280 nm)
<b>Reconstitution</b>	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.
<b>Source</b>	Produced in <i>E. coli</i> . Non-glycosylated, disulfide linked homodimer, containing two identical 113 amino acid chains.

Last modified: May 29, 2024



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)