

# Follistatin (human), (recombinant)

Follistatin is an autocrine acting protein that is expressed by many tissues, but at notably higher levels in the ovary and skin. Follistatin functions to negatively regulate the signaling of a wide variety of TGF- $\beta$  family members (activin, BMPs, myostatin, GDF-11 and TGF- $\beta$ 1). Mechanistically, follistatin works as an antagonist by complexing with TGF- $\beta$  family members to prevent them from interacting with their signaling receptors.

## Ordering Information

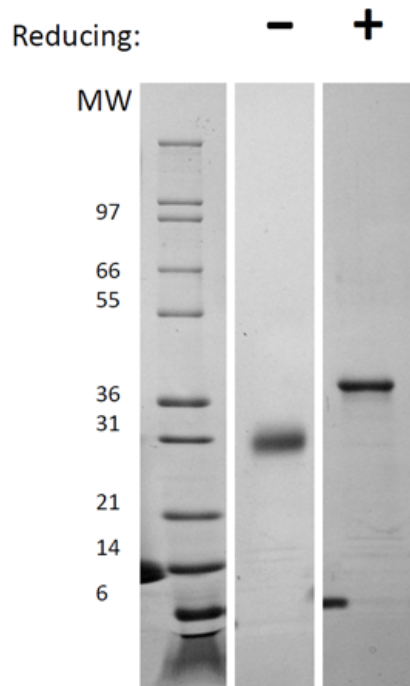
[Order Online »](#)

ALX-201-813-0020	20 $\mu$ g
ALX-201-813-0100	100 $\mu$ g

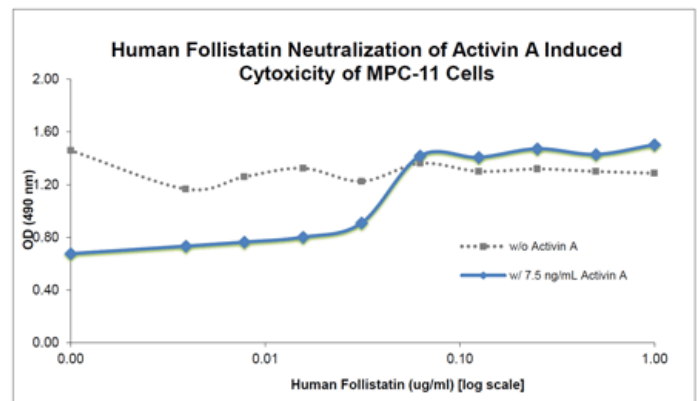
## Manuals, SDS & CofA

[View Online »](#)

- Carrier-free



**Figure 2:** 1µg in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human Follistatin has a predicted MW of 31.7 kDa.



**Figure 1: Human Follistatin Bioactivity Data.** MPC-11 cells were cultured with 0 to 1µg/ml human Follistatin with or without 7.5ng/ml human Activin A. Cell viability was measured after 66 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of human Follistatin was 30-50ng/ml. Complete neutralization is seen by 250ng/ml.

## 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>	FS, Activin-binding protein, FSH-suppressing protein
<b>Appearance</b>	White lyophilized (freeze-dried) powder.
<b>Biological Activity</b>	The activity is determined by the dose-dependent proliferation of mouse BALB/c 3T3 cells which is typically observed at concentrations <0.1ng/ml.
<b>Endotoxin Content</b>	≤1 EU/μg protein measured by kinetic LAL analysis.
<b>Formulation</b>	Lyophilized from 10mM sodium phosphate and 50mM sodium chloride, pH 7.5, sterile filtered.
<b>MW</b>	~31.7kDa
<b>Purity</b>	≥95% (SDS-PAGE using Coomassie Staining)
<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 water 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 protein, containing 289 amino acids.

**UniProt ID** P19883



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)