

Follistatin (mouse), (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- β family members (activin, BMPs, myostatin, GDF-11 and TGF- β 1). Mechanistically, follistatin works as an antagonist by complexing with TGF- β family members to prevent them from interacting with their signaling receptors.

Ordering Information

[Order Online »](#)

ALX-201-814-0020	20 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

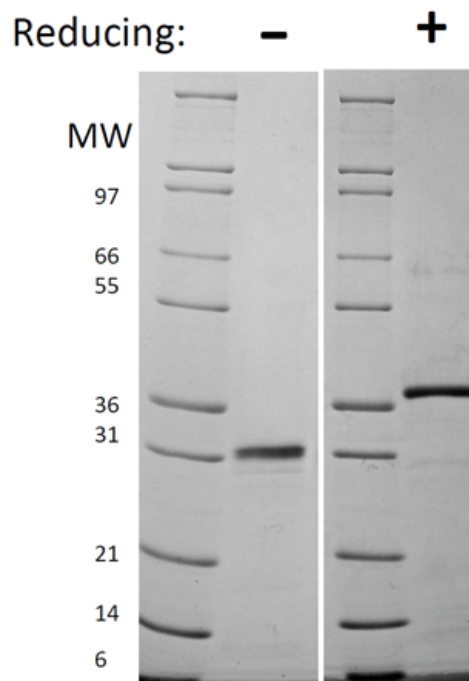


Figure 2: 1 μ g in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Mouse Follistatin has a predicted MW of 31.6 kDa.

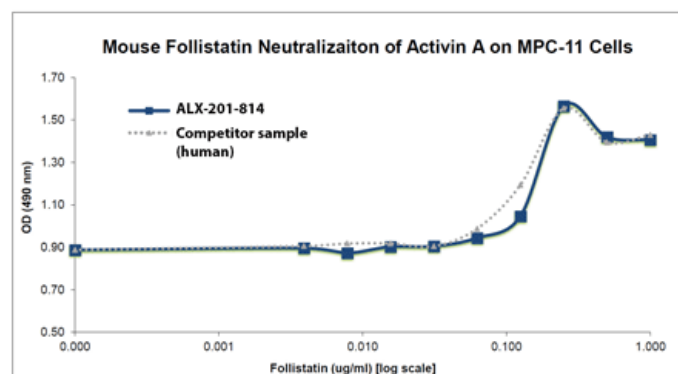


Figure 1: Mouse Follistatin Bioactivity Data. Serial dilutions of mouse Follistatin (starting at 1 μ g/ml) were mixed with 7.5ng/ml human Activin A, then added to MCP-11 cells. After 69 hours cell proliferation was measured and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of mouse Follistatin was between 0.13-0.19 μ g/ml. This value is comparable to the typically expected to be less than 0.3 μ g/ml.

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	FS, Activin-binding protein, FSH-suppressing protein
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the dose-dependent neutralization of 7.5ng/ml human Activin A. Complete neutralization is typically reached at <0.3µg/ml.
Endotoxin Content	≤5 EU/µg protein measured by kinetic LAL analysis.
Formulation	Lyophilized from 10mM Na ₂ PO ₄ , pH 7.5, containing 50mM sodium chloride. Sterile filtered.
MW	~31.6kDa
Purity	≥90% (Reducing and Non-reducing SDS-PAGE)
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 water 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 protein, containing 289 amino acids.

UniProt ID P47931



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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