

Activin-A (human), (recombinant)

Activin A is a member of the TGF- β family of proteins produced by many cell types throughout development. Activins interact with Type I and Type II serine/threonine kinases to signal to SMAD proteins to regulate a variety of functions, including cell proliferation, differentiation, wound healing, apoptosis, metabolism, etc. Activin A is a homodimer of two beta A chains and is not biologically active until the N terminal propeptide is cleaved from each. Human Activin A has 100% amino acid sequence identity to mouse, rat, pig, bovine and cat proteins.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-809-0010	10 μ g
ALX-201-809-0100	100 μ g

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

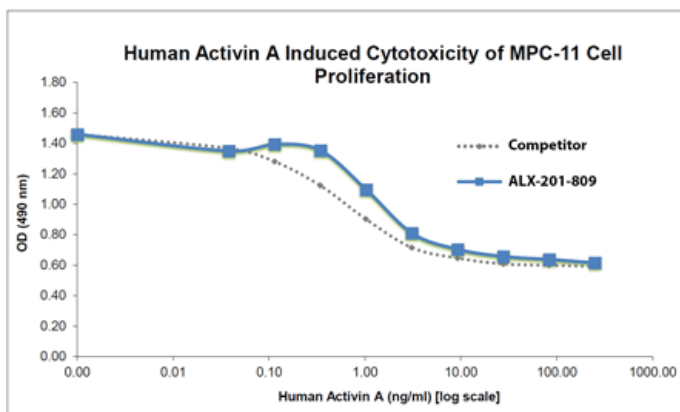


Figure 1: Human Activin A Bioactivity Data. MPC-11 cells were cultured with 0 to 250ng/ml human Activin A. Cell viability was measured after 66 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of human Activin A was 1-1.5ng/ml. This value is comparable to the typical expected range of 1-3ng/ml.

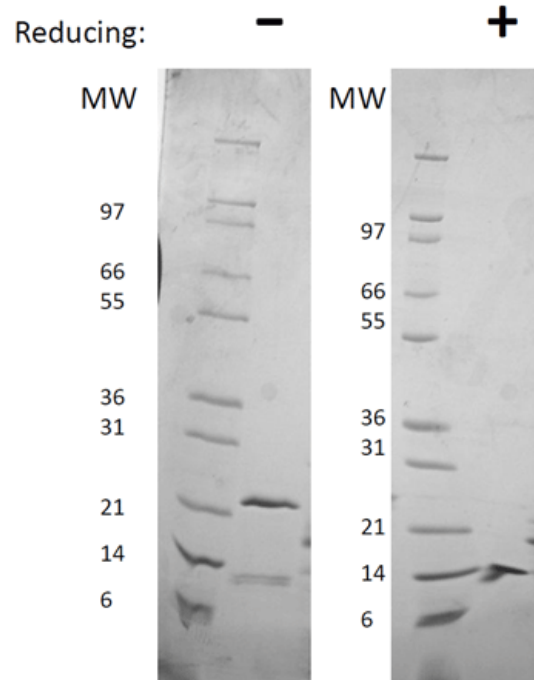


Figure 2: 1µg in each lane (-) non-reducing conditions and (+) reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Human Activin A is a homodimer with a predicted MW of 26.2 kDa.

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	Inhibin β -1, FRP, FSH-releasing protein
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the ability to induce cytotoxicity of MPC-11 cells which is typically observed at concentrations ≤ 10 ng/ml.
Endotoxin Content	≤ 1 EU/ μ g protein measured by kinetic LAL analysis.
Formulation	Lyophilized from a sterile filtered solution containing 0.1% TFA.
MW	~26.2kDa
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 water, which can then be further diluted to other aqueous solutions.
Source	Produced in <i>E. coli</i> . Non-glycosylated homodimer, containing two 117 amino acid chains.
UniProt ID	P08476



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)