# Activin-A (human), (recombinant)

Activin A is a member of the TGF- $\beta$  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**

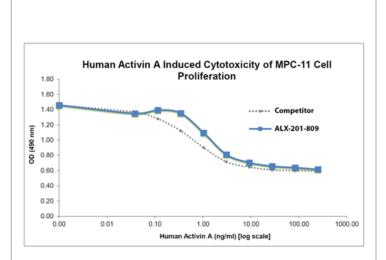
**Order Online** »

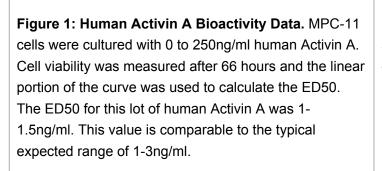
ALX-201-809-0010	10µg
ALX-201-809-0100	100µg

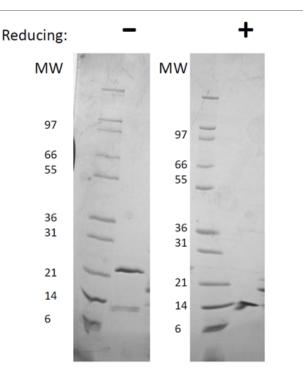
Manuals, SDS & CofA

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 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 cytoxicity

of MPC-11 cells which is typically observed at

concentrations ≤10ng/ml.

Endotoxin Content ≤1 EU/µg protein measured by kinetic LAL analysis.

Formulation Lyophilized from a sterile filtered solution containing 0.1%

TFA.

**MW** ∼26.2kDa

**Purity** ≥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 *E. coli*. Non-glycosylated homodimer,

containing two 117 amino acid chains.

UniProt ID P08476

Last modified: May 29, 2024



info-

eu@enzolifesciences.com