

Epidermal growth factor (rat), (recombinant)

Epidermal Growth Factor (EGF) is a growth factor that stimulates the proliferation of epithelial and epidermal cells. EGF family members are characterized by three intramolecular disulfide bonds and can bind to four different receptor tyrosine kinases known as EGFR/ErbB1, ErbB2, ErbB3, and ErbB4.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-812-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Carrier-free

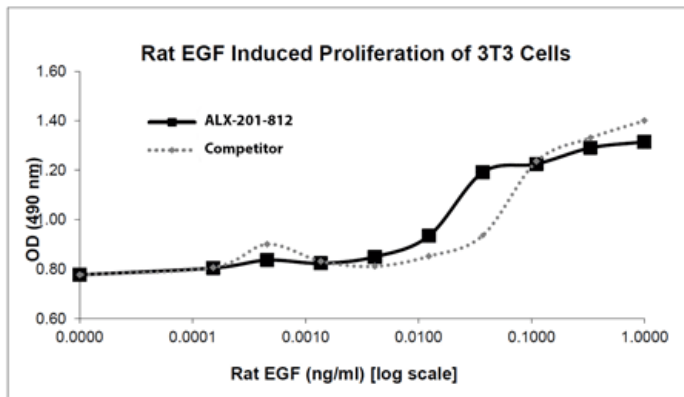


Figure 1: Rat EGF Bioactivity Data. 3T3 cells were cultured with 0 to 1ng/ml Rat EGF. Cell proliferation was measured after 42 hours and the linear portion of the curve was used to calculate the ED50. The ED50 for this lot of Rat EGF was 50-80pg/ml. This value is comparable to the typical expected range of 20-100pg/ml.

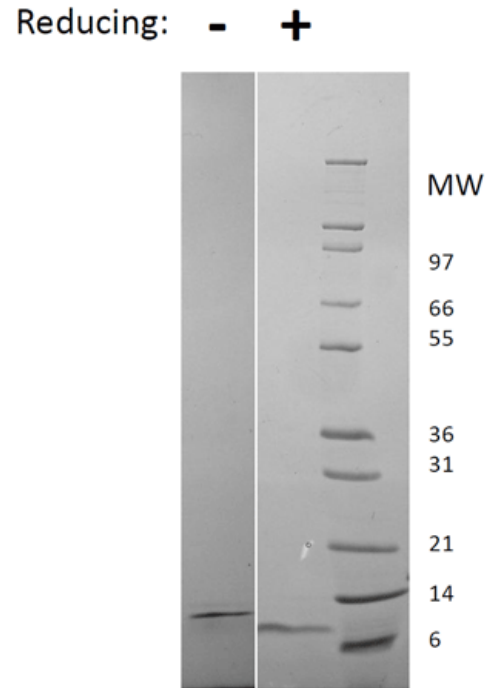


Figure 2: 1µg of protein run under (+) reducing and (-) non-reducing conditions in a 4-20% TRIS-Glycine gel, stained with Coomassie Blue. Rat EGF is predicted to have a MW of 6.3 kDa.

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	EGF, Urogastrone, URG
Appearance	White lyophilized (freeze-dried) powder.
Biological Activity	The activity is determined by the dose-dependent proliferation of mouse BALB/c 3T3 cells which is typically observed at concentrations <0.1ng/ml.
Endotoxin Content	≤1 EU/μg protein measured by kinetic LAL analysis.
Formulation	Lyophilized from 10mM Na ₂ PO ₄ , pH 7.5. Sterile filtered.
MW	~6.3kDa
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 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 54 amino acids, with a molecular weight of 6.3 kDa.
UniProt ID	P07522



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)