

Epidermal growth factor (mouse), (recombinant)

Epidermal Growth Factor (EGF) is a growth factor that stimulates the proliferation of epithelial and epidermal cells. EGF family members can bind to different receptor tyrosine kinases.

Manuals, SDS & CofA

[View Online »](#)

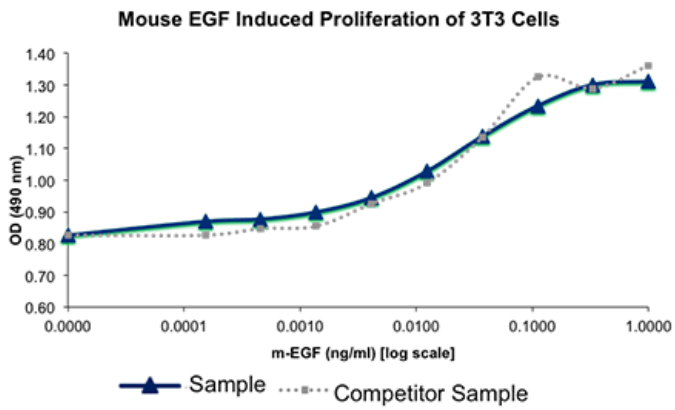


Figure 1: 3T3 cells were cultured with 0 to 1 ng/mL Mouse EGF. Cell proliferation was measured after 44 hours and the linear portion of the curve can be used to calculate the ED50.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Urogastrone, URG, EGF
Endotoxin Content	≤ 0.05 EU/μg protein (LAL test)
Formulation	Lyophilized from 0.2 μm filtered solution containing 10mM sodium phosphate, pH 7.5
MW	~6 kDa
Purity	≥95% (Coomassie staining)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . A non-glycosylated protein containing 53 amino acids.
UniProt ID	P01132

Last modified: May 29, 2024



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