

Epidermal growth factor (human), (recombinant)

Recombinant human EGF that regulates cell growth, proliferation, and differentiation.

Human epidermal growth factor (EGF, also known as beta-urogastrone) is a 53 amino acid peptide that regulates cell growth, proliferation, and differentiation. EGF is an integral membrane protein that is cleaved and released into the extracellular space where it acts as a hormone. EGF binds cell surface EGF receptors (EGFR), initiating signaling cascades through activation of EGFR tyrosine kinase activity.

Ordering Information

[Order Online »](#)

ADI-908-060-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name EGF

Application Notes Western blot control.

Formulation EGF was lyophilized from a concentrated (1mg/mL) solution containing PBS pH-7.4.

Purity ≥95% (SDS-PAGE)

Source Produced in *E. coli*.

UniProt ID P01133

Last modified: October 9, 2025



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)