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



eu@enzolifesciences.com