

Recombinant Human Epidermal Growth Factor (EGF)

Cat. No.: 13247-051

Size: 100 µg

Store at -20°C.

Description:

Recombinant Human Epidermal Growth Factor, from *E.coli*, is a 6.2 kDa protein containing 53 amino-acid residues. It is a polypeptide growth factor which stimulates the proliferation of a wide range of epidermal and epithelial cells (1-4). For most *in vitro* applications, EGF exerts its biological activity in the concentration range of 0.5 to 25 ng/ml. Recombinant Human EGF is provided as a lyophilized protein obtained from a membrane-filtered (0.22 µm) solution of distilled water with no additives.

Directions for Use:

1. Reconstitute in distilled water, aqueous buffers or 10 mM acetic acid to a concentration of 1 µg/µl. **Note: Do not store in glass;** store in polypropylene vial. Store at -20°C reconstituted for up to six months. Avoid repeated freeze-thaw cycles.
2. The reconstituted solution can be further diluted in distilled water or aqueous buffers to working concentration.

Note: To avoid adsorptive loss, dilute solutions should be supplemented with Bovine Serum Albumin (BSA) at a concentration of not less than 0.1%.

Doc. Rev.: 05/24/01

Quality Control:

- Purity: $\geq 98\%$ (by SDS-PAGE)
- Endotoxin: ≤ 0.1 ng/ μ g of EGF
- Biological Activity: $ED_{50} \leq 2$ ng/ml (determined by dose-dependent stimulation of thymidine uptake by BALB/c 3T3)
- Specific Activity: $\geq 5 \times 10^5$ units/mg

References:

1. Cohen, S. (1960) *Proc. Natl. Acad. Sci.* 46, 302.
2. Levi-Montalcini, R. *et al.* (1960) *Ann. Rev. Acad. Sci.* 85, 324.
3. Savage, C.R. (1973) *J. Biol. Chem.* 247, 7612.
4. Messing, E.M. *et al.* (1987) *Cancer, Res.* 47, 2230.