

---

## EGF Polyclonal Antibody

(Catalog # A70851)

---

### Background

This gene encodes a member of the epidermal growth factor superfamily. The encoded preproprotein is proteolytically processed to generate the 53-amino acid epidermal growth factor peptide. This protein acts a potent mitogenic factor that plays an important role in the growth, proliferation and differentiation of numerous cell types. This protein acts by binding with high affinity to the cell surface receptor, epidermal growth factor receptor. Defects in this gene are the cause of hypomagnesemia type 4. Dysregulation of this gene has been associated with the growth and progression of certain cancers. Alternative splicing results in multiple transcript variants, at least one of which encodes a preproprotein that is proteolytically processed.

### Description

EGF Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

P01133

### Purification

Affinity Purification

### Immunogen

Recombinant protein of human EGF

### Storage

Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

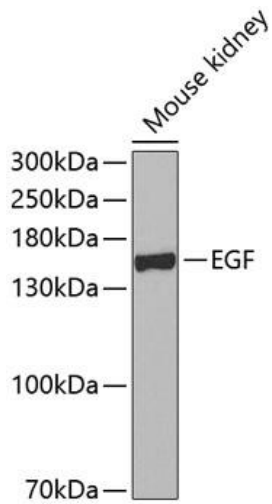
### Alternative Names

EGF; HOMG4; URG; epidermal growth factor

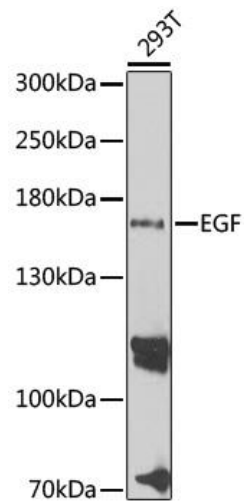
### Application

DB, WB, Recommended dilution: DB 1:500 - 1:2000, WB 1:500 - 1:2000

---



Western blot analysis of extracts of mouse kidney, using EGF Polyclonal Antibody at 1:500 dilution.  
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.  
Lysates/proteins: 25ug per lane.  
Blocking buffer: 3% nonfat dry milk in TBST.  
Exposure time: 90s.



Western blot analysis of extracts of 293T cells, using EGF Polyclonal Antibody at 1:1000 dilution.  
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.  
Lysates/proteins: 25ug per lane.  
Blocking buffer: 3% nonfat dry milk in TBST.  
Exposure time: 10s.