

---

## LIFR Polyclonal Antibody

(Catalog # A69575)

---

### Background

Signal-transducing molecule. May have a common pathway with IL6ST. The soluble form inhibits the biological activity of LIF by blocking its binding to receptors on target cells.

### Description

LIFR Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, pH 7.4

### Specificity

Human, Rat

### Isotype

IgG

### Uniprot ID

P42702

### Purification

Protein G purified

### Immunogen

Recombinant Human Leukemia inhibitory factor receptor protein (915-1086AA)

### Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

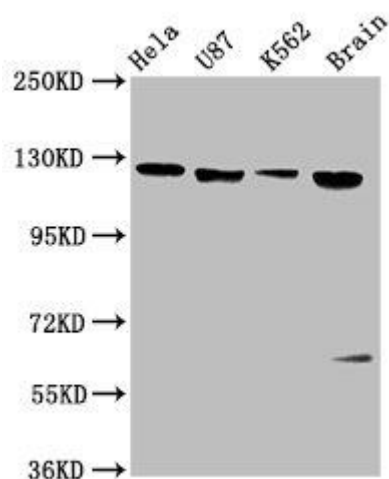
### Alternative Names

Leukemia inhibitory factor receptor, LIF receptor, LIF-R, CD118, LIFR

### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:500-1:1000, IF:1:200-1:500

---



#### Western Blot

Positive WB detected in: HeLa whole cell lysate, U87 whole cell lysate, K562 whole cell lysate, Rat brain tissue

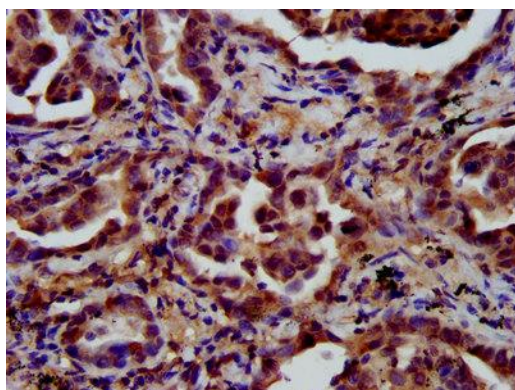
All lanes: LIFR antibody at 3.4ug/ml

#### Secondary

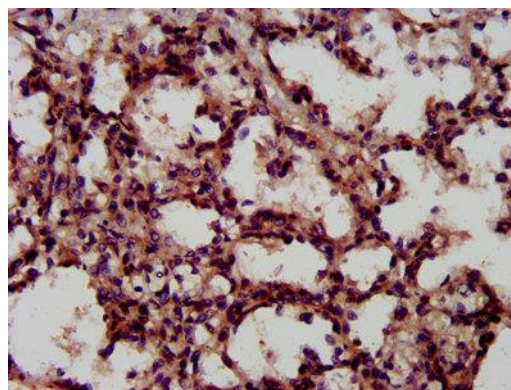
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 124 KDa

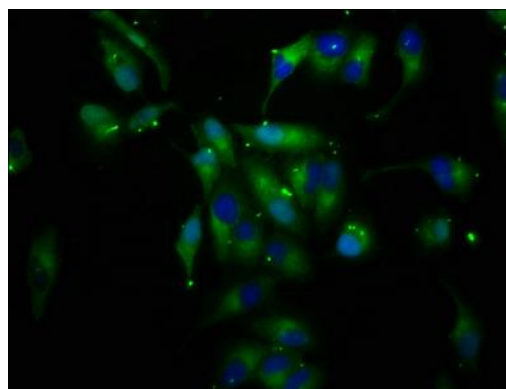
Observed band size: 124 KDa



IHC image of LIFR Polyclonal Antibody diluted at 1:600 and staining in paraffin-embedded human lung cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of LIFR Polyclonal Antibody diluted at 1:600 and staining in paraffin-embedded human lung tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HeLa cells with LIFR Polyclonal Antibody at 1:200, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).