

HBEGF Polyclonal Antibody

(Catalog # A73040)

Description

HBEGF Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q99075

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 20-160 of human HBEGF (NP_001936.1).

Storage

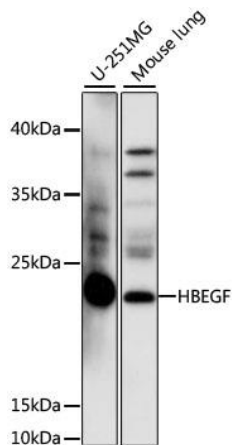
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

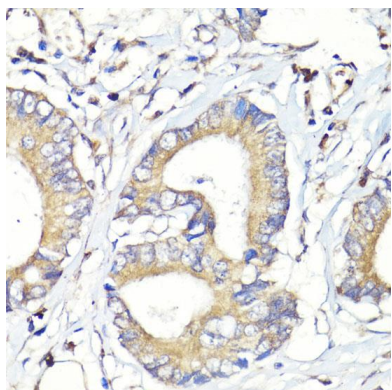
HBEGF; DTR; DTS; DTSF; HEGFL; heparin binding EGF like growth factor

Application

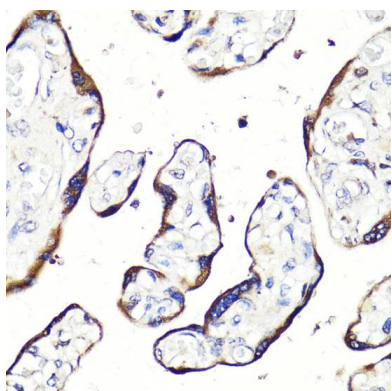
WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IF 1:50 - 1:200



Western blot analysis of extracts of various cell lines, using HBEGF 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: 90s.



Immunohistochemistry of paraffin-embedded human colon carcinoma using HBEGF Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human placenta using HBEGF Polyclonal Antibody at dilution of 1:100 (40x lens).