
IGF1 Polyclonal Antibody

(Catalog #A70939)

Background

The protein encoded by this gene is similar to insulin in function and structure and is a member of a family of proteins involved in mediating growth and development. The encoded protein is processed from a precursor, bound by a specific receptor, and secreted. Defects in this gene are a cause of insulin-like growth factor I deficiency. Alternative splicing results in multiple transcript variants encoding different isoforms that may undergo similar processing to generate mature protein.

Description

IGF1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.09% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P05019

Purification

Affinity Purification

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 50-150 of human IGF1 (NP_000609.1).

Storage

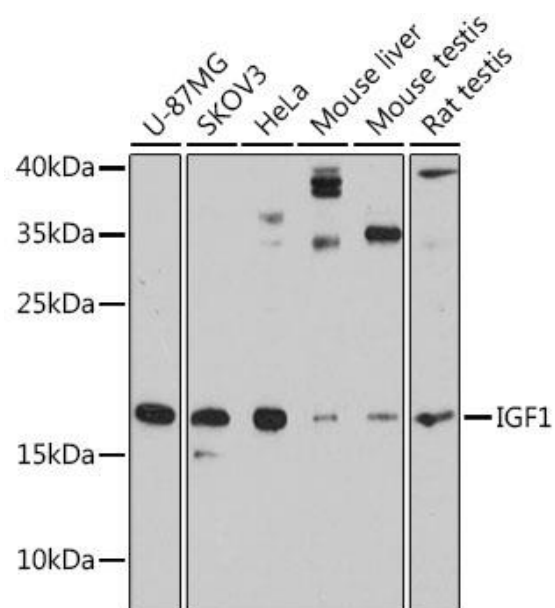
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

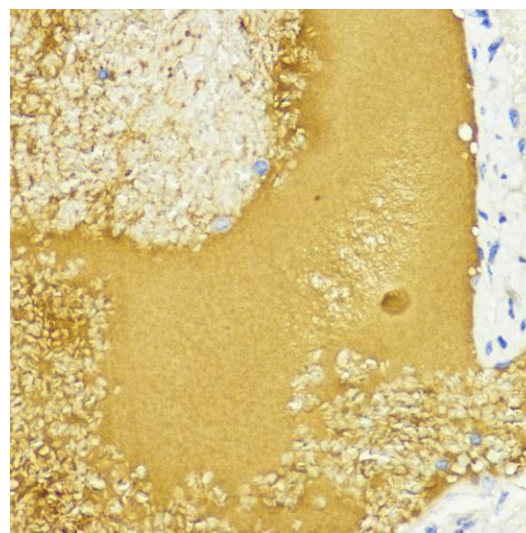
IGF1; IGF-I; IGF1; MGF; insulin-like growth factor I

Applications

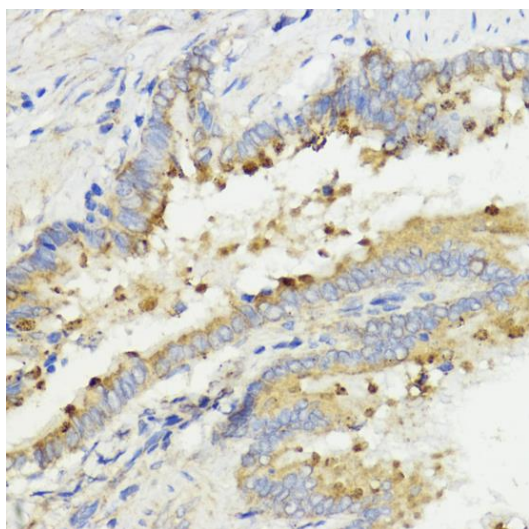
WB, IHC, ELISA; Recommended dilution: WB 1:1500 - 1:3000, IHC 1:50 - 1:100, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



Western blot analysis of extracts of various cell lines, using IGF1 Polyclonal Antibody at 1:3000 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 mouse liver using IGF1 Polyclonal Antibody at dilution of 1:100 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.



Immunohistochemistry of paraffin-embedded rat lung using IGF1 Polyclonal Antibody at dilution of 1:100 (40x lens). Microwave antigen retrieval performed with 0.01M PBS Buffer (pH 7.2) prior to IHC staining.