

IFIT3 Polyclonal Antibody

(Catalog #A71627)

Description

IFIT3 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

O14879

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 128-331 of mouse IFIT3 (NP_034631.1).

Storage

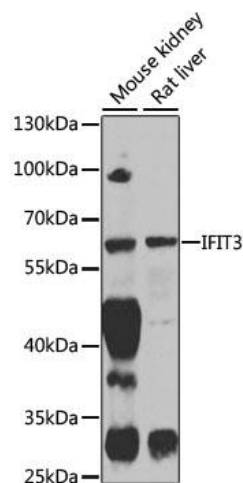
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

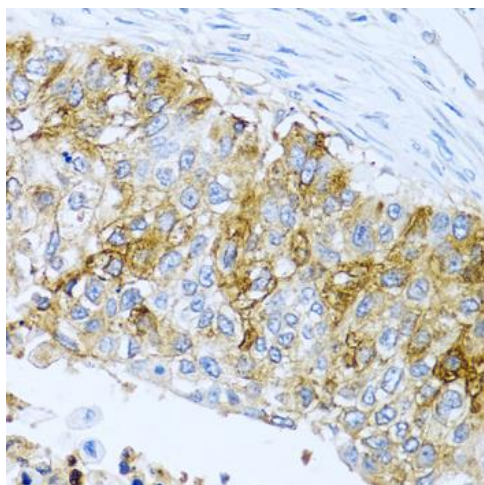
IFIT3; CIG-49; GARG-49; IFI60; IFIT4; IRG2; ISG60; P60; RIG-G; cig41; interferon-induced protein with tetratricopeptide repeats 3

Application

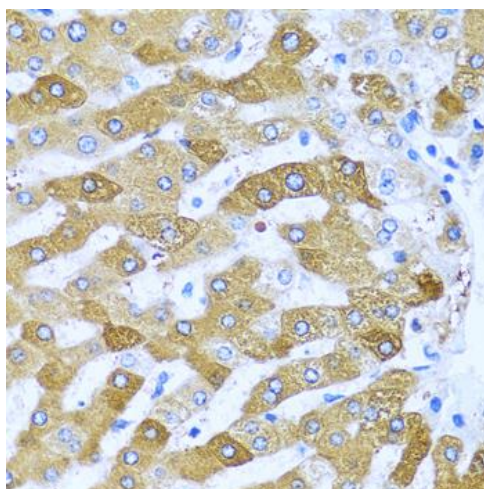
ELISA, WB, IHC; Recommended dilution: WB 1:500 – 1:1000, IHC 1:100 – 1:200



Western blot analysis of extracts of various cell lines, using IFIT3 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: 60s.



Immunohistochemistry of paraffin-embedded human lung cancer using IFIT3 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human liver using IFIT3 Polyclonal Antibody at dilution of 1:100 (40x lens).