
IFIT5 Polyclonal Antibody

(Catalog A59462)

Background

Interferon-induced RNA-binding protein that specifically binds single-stranded RNA bearing a 5'-triphosphate group (PPP-RNA), thereby acting as a sensor of viral single-stranded RNAs. Single-stranded PPP-RNAs, which lack 2'-O-methylation of the 5' cap and bear a 5'-triphosphate group instead, are specific from viruses, providing a molecular signature to distinguish between self and non-self mRNAs by the host during viral infection. Directly binds PPP-RNA in a non-sequence-specific manner. Also recognizes and binds tRNAs.

Description

IFIT5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q13325

Purification

Protein G purified

Immunogen

Recombinant Human Interferon-induced protein with tetratricopeptide repeats 5 protein (107-320AA)

Storage

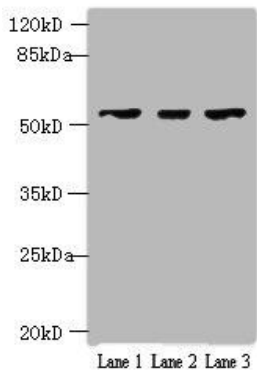
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

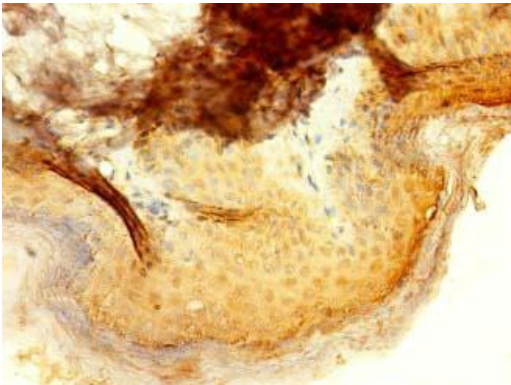
Interferon-induced protein with tetratricopeptide repeats 5, IFIT-5, IFIT5, ISG58, RI58, Interferon-induced 58 kDa protein, Retinoic acid- and interferon-inducible 58 kDa protein, P58

Application

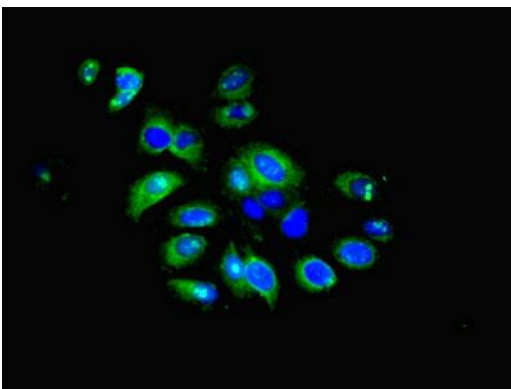
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western blot
All lanes: IFIT5 Polyclonal Antibody at 2ug/ml
Lane 1: A375 whole cell lysate
Lane 2: Jurkat whole cell lysate
Lane 3: HepG2 whole cell lysate
Secondary
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 56, 51 kDa
Observed band size: 56 kDa



Immunohistochemistry of paraffin-embedded human skin tissue using IFIT5 Polyclonal Antibody at dilution 1:100



Immunofluorescent analysis of HepG2 cells using IFIT5 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)