
Ifne Polyclonal Antibody

(Catalog # A55174)

Background

Type I interferon required for maintaining basal levels of IFN-regulated genes, including 2'-5'-oligoadenylate synthetase, IRF7 and ISG15, in the female reproductive tract. Directly mediates protection against viral, including HSV-2, and bacterial, including Chlamydia muridarum, genital infections.

Description

Ifne Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

Q80ZF2

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse Interferon epsilon protein (22-192AA)

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

IFN-epsilon, Interferon epsilon-1, Interferon tau-1, Ifne, Ifne1, Ifnt1

Application

ELISA
