
ERVK-6 Polyclonal Antibody

(Catalog A65250)

Background

Retroviral envelope proteins mediate receptor recognition and membrane fusion during early infection. Endogenous envelope proteins may have kept, lost or modified their original function during evolution. This endogenous envelope protein has lost its original fusogenic properties.

Description

ERVK-6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q69384

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Endogenous retrovirus group K member 6 Env polyprotein (90-632AA)

Storage

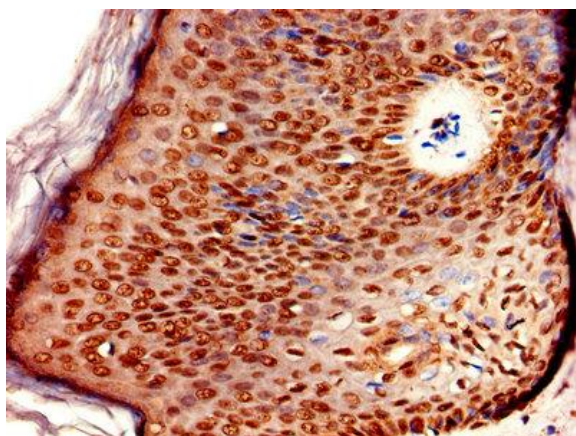
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

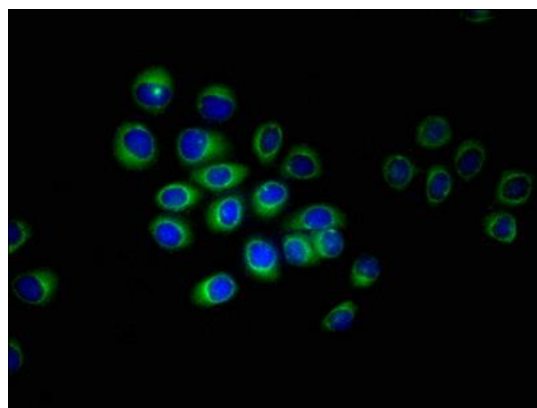
Endogenous retrovirus group K member 6 Env polyprotein, EnvK2 protein, Envelope polyprotein, HERV-K(C7) envelope protein, HERV-K(HML-2.HOM) envelope protein, HERV-K108 envelope protein, HERV-K_7p22.1 provirus ancestral Env polyprotein, ERVK-6, ERVK6

Application

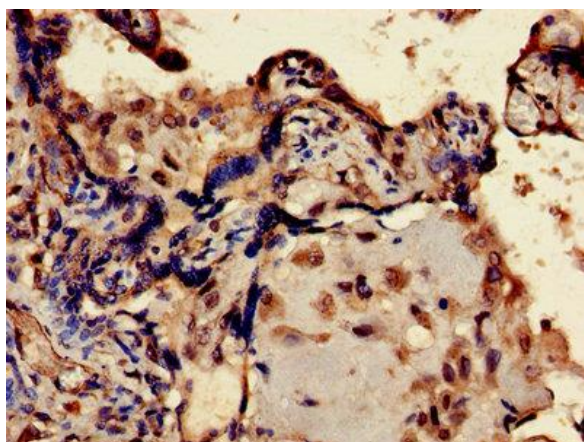
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human skin tissue using ERVK-6 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of A549 cells using ERVK-6 Polyclonal Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human placenta tissue using ERVK-6 Polyclonal Antibody at dilution of 1:100