

IL24 Polyclonal Antibody

(Catalog # A64634)

Background

Has antiproliferative properties on melanoma cells and may contribute to terminal cell differentiation.

Description

IL24 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

Q13007

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Interleukin-24 protein (52-207AA)

Storage

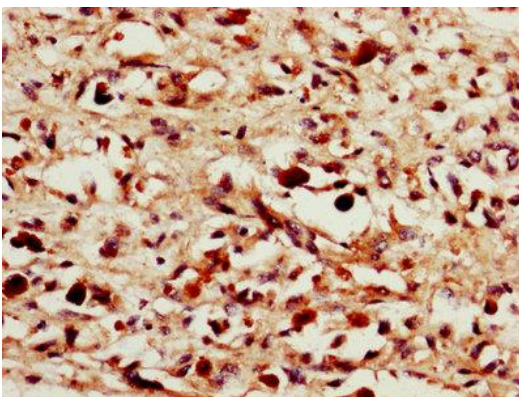
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

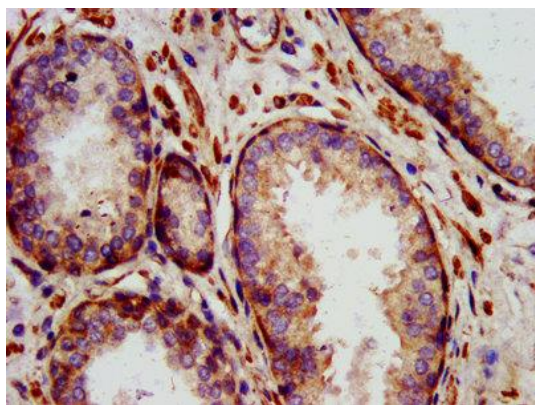
Interleukin-24, IL-24, Melanoma differentiation-associated gene 7 protein, MDA-7, Suppression of tumorigenicity 16 protein, IL24, MDA7, ST16

Application

ELISA, IHC; Recommended dilution: IHC:1:1000-1:2000



IHC image of IL24 Polyclonal Antibody diluted at 1:1400 and staining in paraffin-embedded human melanoma. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of IL24 Polyclonal Antibody diluted at 1:1400 and staining in paraffin-embedded human prostate cancer. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.