
DPY30 Polyclonal Antibody

(Catalog #A58846)

Background

As part of the MLL1/MLL complex, involved in the methylation of histone H3 at 'Lys-4', particularly trimethylation. Histone H3 'Lys-4' methylation represents a specific tag for epigenetic transcriptional activation. May play some role in histone H3 acetylation. In a teratocarcinoma cell, plays a crucial role in retinoic acid-induced differentiation along the neural lineage, regulating gene induction and H3 'Lys-4' methylation at key developmental loci. May also play an indirect or direct role in endosomal transport.

Description

DPY30 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

Q9C005

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein dpy-30 homolog protein (1-99AA)

Storage

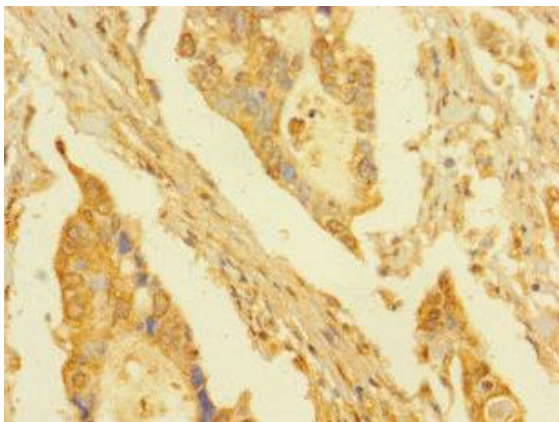
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

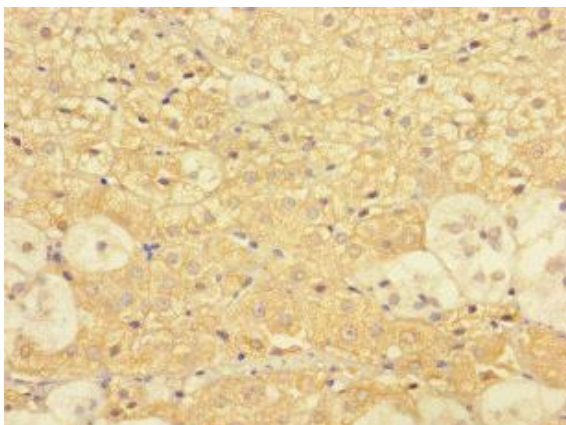
Protein dpy-30 homolog, DPY30, Dpy-30-like protein, Dpy-30L

Application

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



Immunohistochemistry of paraffin-embedded human pancreatic cancer using DPY30 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using DPY30 Polyclonal Antibody at dilution of 1:100