

ALKBH3 Polyclonal Antibody

(Catalog # A57365)

Background

Dioxygenase that repairs alkylated DNA containing 1-methyladenine (1meA) and 3-methylcytosine (3meC) by oxidative demethylation. Has a strong preference for single-stranded DNA. Able to process alkylated 3mC within double-stranded regions via its interaction with ASCC3, which promotes DNA unwinding to generate single-stranded substrate needed for ALKBH3. May also act on RNA. Requires molecular oxygen, alpha-ketoglutarate and iron.

Description

ALKBH3 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

Q96Q83

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Alpha-ketoglutarate-dependent dioxygenase alkB homolog 3 protein (1-170AA)

Storage

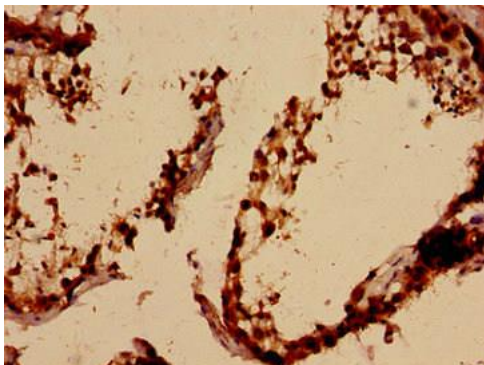
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

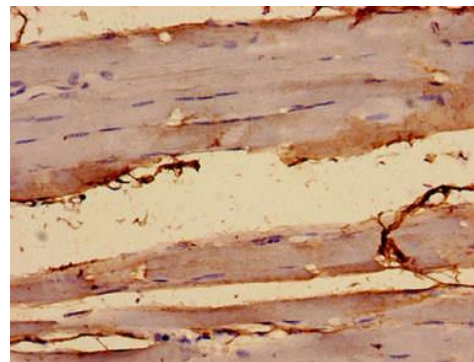
Alkylated DNA repair protein alkB homolog 3, DEPC-1, Prostate cancer antigen 1, ABH3, DEPC1, ALKBH3

Application

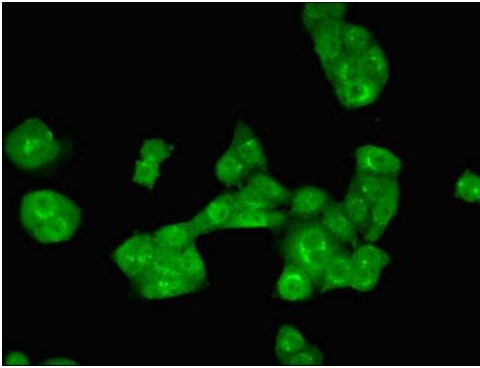
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry analysis of human testis tissue using ALKBH3 Polyclonal Antibody at dilution of 1:100



Immunocytochemistry analysis of human skeletal muscle tissue using ALKBH3 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of PC3 cells using ALKBH3 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)