
NANOS2 Polyclonal Antibody

(Catalog #A63006)

Background

Plays a key role in the sexual differentiation of germ cells by promoting the male fate but suppressing the female fate. Represses the female fate pathways by suppressing meiosis, which in turn results in the promotion of the male fate. Maintains the suppression of meiosis by preventing STRA8 expression, which is required for premeiotic DNA replication, after CYP26B1 is decreased. Regulates the localization of the CCR4-NOT deadenylation complex to P-bodies and plays a role in recruiting the complex to trigger the degradation of mRNAs involved in meiosis. Required for the maintenance of the spermatogonial stem cell population. Not essential for the assembly of P-bodies but is required for the maintenance of their normal state.

Description

NANOS2 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

P60321

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Nanos homolog 2 protein (1-138AA)

Storage

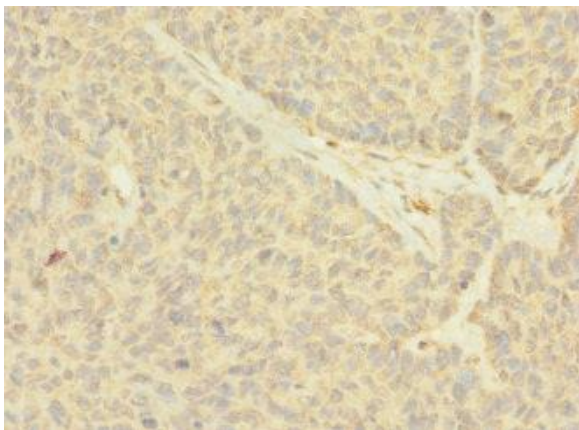
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

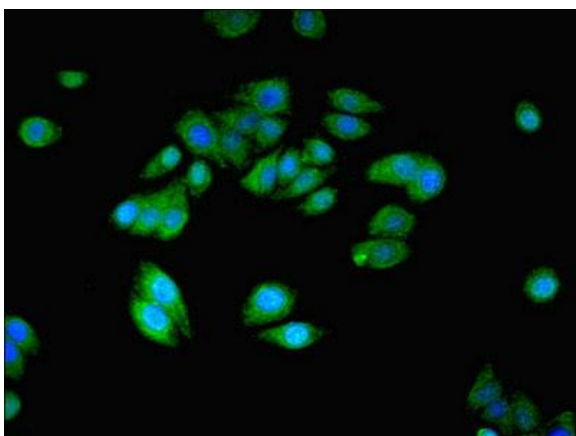
Nanos homolog 2, NOS-2, NANOS2, NOS2

Application

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



Immunohistochemistry of paraffin-embedded human ovarian cancer using NANOS2 Polyclonal Antibody at dilution 1: 100



Immunofluorescent analysis of HepG2 cells using NANOS2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)