
ZNF281 Polyclonal Antibody

(Catalog #A63844)

Background

Transcription repressor that plays a role in regulation of embryonic stem cells (ESCs) differentiation. Required for ESCs differentiation and acts by mediating autorepression of NANOG in ESCs: binds to the NANOG promoter and promotes association of NANOG protein to its own promoter and recruits the NuRD complex, which deacetylates histones. Not required for establishment and maintenance of ESCs. Represses the transcription of a number of genes including GAST, ODC1 and VIM. Binds to the G-rich box in the enhancer region of these genes.

Description

ZNF281 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

Q9Y2X9

Purification

Antigen Affinity purified

Immunogen

Recombinant Human Zinc finger protein 281 protein (1-300AA)

Storage

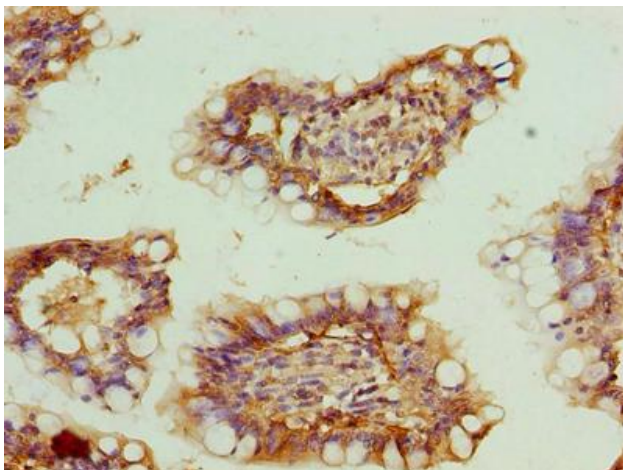
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Zinc finger protein 281, GC-box-binding zinc finger protein 1, Transcription factor ZBP-99, Zinc finger DNA-binding protein 99, ZNF281, GZP1, ZBP99

Application

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



Immunohistochemistry of paraffin-embedded human small intestine tissue using ZNF281 Polyclonal Antibody at dilution of 1:100