
CXXC5 Polyclonal Antibody

(Catalog #A65418)

Background

May indirectly participate in activation of the NF-kappa-B and MAPK pathways. Acts as a mediator of BMP4-mediated modulation of canonical Wnt signaling activity in neural stem cells. Required for DNA damage-induced ATM phosphorylation, p53 activation and cell cycle arrest. Involved in myelopoiesis. Transcription factor. Binds to the oxygen responsive element of COX4I2 and represses its transcription under hypoxia conditions (4% oxygen), as well as normoxia conditions (20% oxygen) (PubMed:23303788). May repress COX4I2 transactivation induced by CHCHD2 and RBPJ (PubMed:23303788).

Description

CXXC5 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

Q7LFL8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human CXXC-type zinc finger protein 5 protein (1-227AA)

Storage

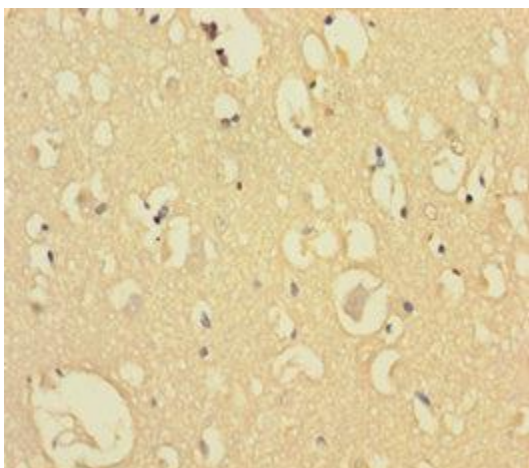
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

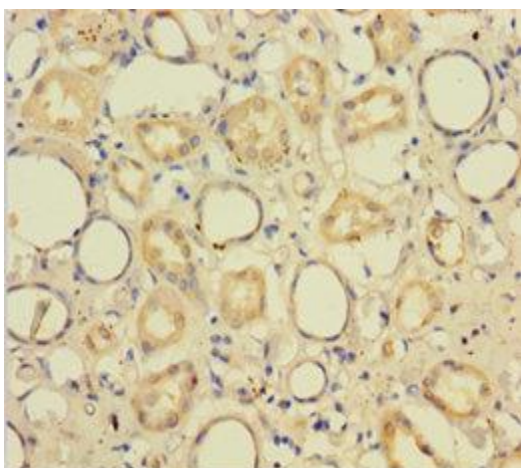
CXXC-type zinc finger protein 5, CF5, Putative MAPK-activating protein PM08, Putative NF-kappa-B-activating protein 102, Retinoid-inducible nuclear factor, RINF, CXXC5, HSPC195, TCCCIA00297

Application

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



Immunohistochemistry of paraffin-embedded human brain tissue using CXXC5 Polyclonal Antibody at dilution 1:100



Immunohistochemistry of paraffin-embedded human kidney tissue using CXXC5 Polyclonal Antibody at dilution 1:100