

HOPX Polyclonal Antibody

(Catalog # A59410)

Background

Atypical homeodomain protein which does not bind DNA and is required to modulate cardiac growth and development. Acts via its interaction with SRF, thereby modulating the expression of SRF-dependent cardiac-specific genes and cardiac development. Prevents SRF-dependent transcription either by inhibiting SRF binding to DNA or by recruiting histone deacetylase (HDAC) proteins that prevent transcription by SRF. Overexpression causes cardiac hypertrophy. May act as a tumor suppressor.

Description

HOPX 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

Q9BPY8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Homeodomain-only protein (1-73AA)

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

Homeodomain-only protein, HOPX, HOD, HOP, LAGY, NECC1, OB1, Lung cancer-associated Y protein, Not expressed in choriocarcinoma protein 1, Odd homeobox protein 1

Application

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



Immunohistochemistry of paraffin-embedded human heart tissue using HOPX Polyclonal Antibody at dilution of 1:100