
ATOH1 Polyclonal Antibody, FITC Conjugated

(Catalog # A56016)

Background

Transcriptional regulator. Activates E box-dependent transcription in collaboration with TCF3/E47, but the activity is completely antagonized by the negative regulator of neurogenesis HES1. Plays a role in the differentiation of subsets of neural cells by activating E box-dependent transcription.

Description

ATOH1 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q92858

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein atonal homolog 1 protein (1-354AA)

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

Class A basic helix-loop-helix protein 14, bHLHa14, Helix-loop-helix protein hATH-1, hATH1, ATH1, BHLHA14, ATOH1

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
