

TLE4 Polyclonal Antibody

(Catalog # A68811)

Background

Transcriptional corepressor that binds to a number of transcription factors. Inhibits the transcriptional activation mediated by PAX5, and by CTNNB1 and TCF family members in Wnt signaling. The effects of full-length TLE family members may be modulated by association with dominant-negative AES. Essential for the transcriptional repressor activity of SIX3 during retina and lens development and for SIX3 transcriptional auto-repression.

Description

TLE4 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

Q04727

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transducin-like enhancer protein 4 protein (284-394AA)

Storage

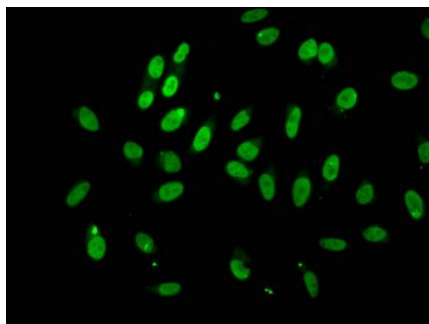
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transducin-like enhancer protein 4, Grg-4, Groucho-related protein 4, TLE4, GRG4, KIAA1261

Application

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



Immunofluorescence staining of HeLa cells with TLE4 Polyclonal Antibody at 1:100, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).