

STOX1 Polyclonal Antibody

(Catalog # A65730)

Background

Involved in regulating the levels of reactive oxidative species and reactive nitrogen species and in mitochondrial homeostasis in the placenta (PubMed:24738702). Required for regulation of inner ear epithelial cell proliferation via the AKT signaling pathway.

Description

STOX1 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

Q6ZVD7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Storkhead-box protein 1 protein (559-858AA)

Storage

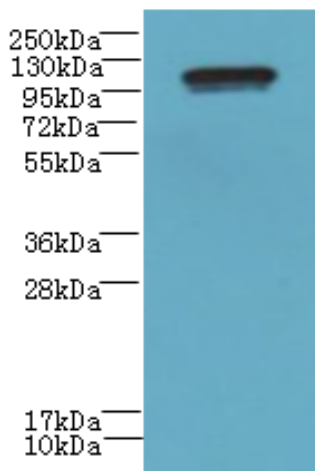
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Storkhead-box protein 1, Winged-helix domain-containing protein, STOX1, C10orf24

Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



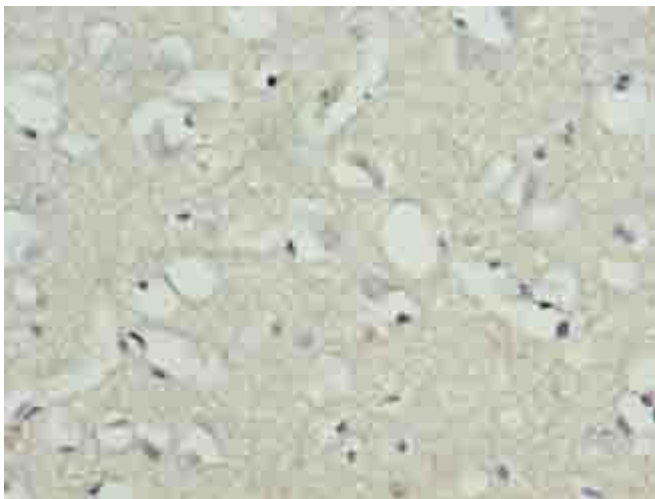
Western blot

All lanes: STOX1 Polyclonal Antibody at 8ug/ml+HepG-2 whole cell lysate

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 111 kDa

Observed band size: 111kDa



Immunohistochemistry of paraffin-embedded human brain tissue using STOX1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human skin tissue using STOX1 Polyclonal Antibody at dilution of 1:100