
DHRS3 Polyclonal Antibody

(Catalog #A62410)

Background

Catalyzes the reduction of all-trans-retinal to all-trans-retinol in the presence of NADPH.

Description

DHRS3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O75911

Purification

Affinity Purified

Immunogen

Recombinant Human Short-chain dehydrogenase/reductase 3 protein (30-169AA)

Storage

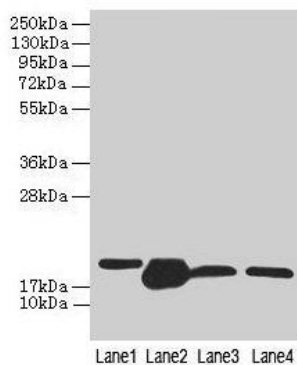
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Short-chain dehydrogenase/reductase 3, DD83.1, Retinal short-chain dehydrogenase/reductase 1, retSDR1, Retinol dehydrogenase 17, Short chain dehydrogenase/reductase family 16C member 1, DHRS3, RDH17, SDR16C1, UNQ2424/PRO4983

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

All lanes: DHRS3 antibody at 0.8ug/ml

Lane 1: A375 whole cell lysate

Lane 2: Mouse liver tissue

Lane 3: Mouse spleen tissue

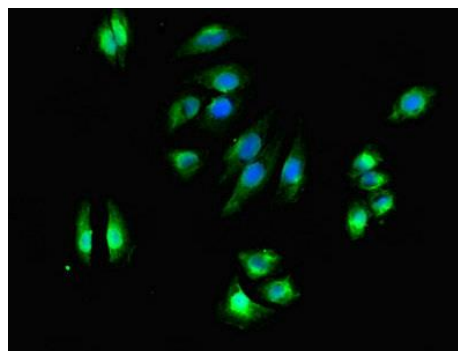
Lane 4: HL60 whole cell lysate

Secondary

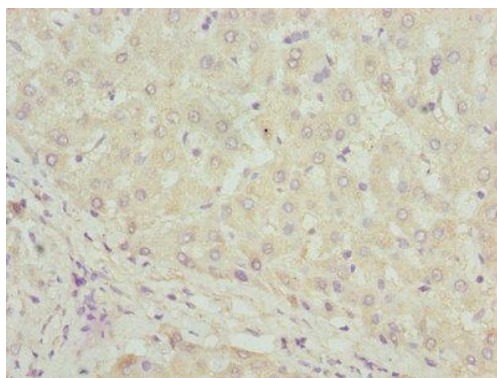
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 34, 19 kDa

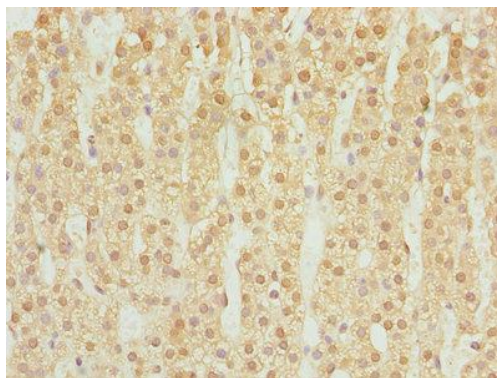
Observed band size: 19 kDa



Immunofluorescent analysis of A549 cells using DHRS3 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human liver tissue using DHRS3 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using DHRS3 Antibody at dilution of 1:100