

RS1 Polyclonal Antibody

(Catalog # A57129)

Background

May be active in cell adhesion processes during retinal development.

Description

RS1 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

O15537

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Retinoschisin protein (24-224AA)

Storage

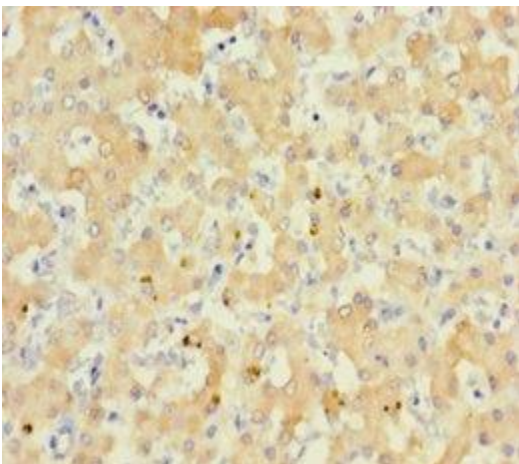
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

X-linked juvenile retinoschisis protein, XLRS1, RS1

Application

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



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