

G0S2 Polyclonal Antibody

(Catalog # A50229)

Background

Promotes apoptosis by binding to BCL2, hence preventing the formation of protective BCL2-BAX heterodimers.

Description

G0S2 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

P27469

Purification

Protein G purified

Immunogen

Recombinant Human G0/G1 switch protein 2 protein (1-103AA)

Storage

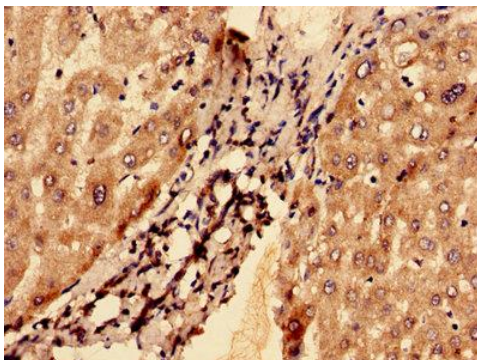
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

G0/G1 switch protein 2, G0/G1 switch regulatory protein 2, Putative lymphocyte G0/G1 switch gene, G0S2

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200



Immunohistochemistry analysis of human liver tissue using G0S2 Antibody at dilution of 1:100