

HBB Polyclonal Antibody

(Catalog # A50281)

Background

Involved in oxygen transport from the lung to the various peripheral tissues. LVV-hemorphin-7 potentiates the activity of bradykinin, causing a decrease in blood pressure. Spinorphin: functions as an endogenous inhibitor of enkephalin-degrading enzymes such as DPP3, and as a selective antagonist of the P2RX3 receptor which is involved in pain signaling, these properties implicate it as a regulator of pain and inflammation.

Description

HBB Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

P68871

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Hemoglobin subunit beta protein (1-147AA)

Storage

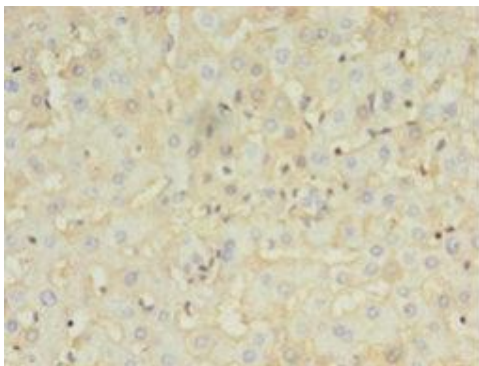
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

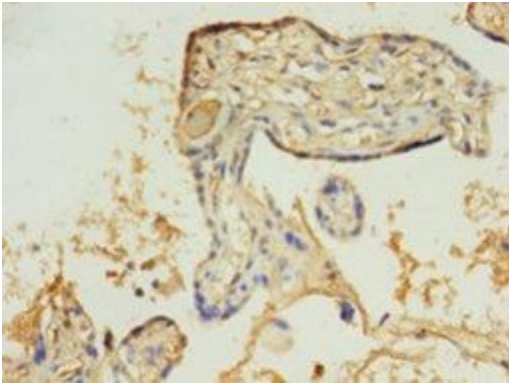
Beta-globin, Hemoglobin beta chain, HBB

Application

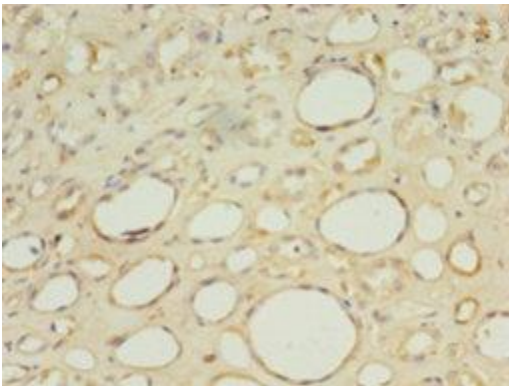
ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:200-1:500



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



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



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