

Mouse Hemoglobin Polyclonal Antibody

(Catalog # A57573)

Background

Hemoglobin is involved in oxygen transport from the lung to the various peripheral tissues. The alpha (HBA) and beta (HBB) loci determine the structure of the 2 types of polypeptide chains in adult Hemoglobin. The normal adult Hemoglobin tetramer consists of two alpha chains and two beta chains. Mutant beta globin causes sickle cell anemia. Absence of beta chain causes beta zero thalassemia. Reduced amounts of detectable beta globin causes beta plus thalassemia.

Description

Mouse Hemoglobin Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Purification

>95%, Protein G purified

Immunogen

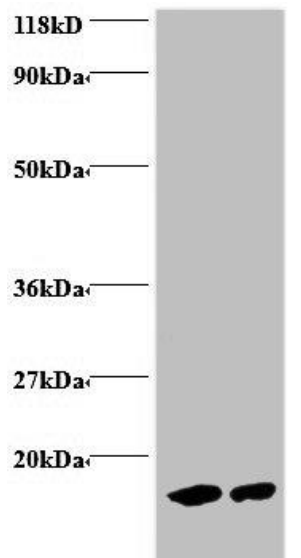
Mouse Hemoglobin protein (Native Protein)

Storage

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

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot
All lanes: Mouse Hemoglobin Polyclonal Antibody at 2ug/ml+
Mouse serum
Lane 1: Mouse serum at 1:100
Lane 2: Mouse serum at 1:1000
Secondary
Goat polyclonal to Rabbit IgG at 1/15000 dilution
Predicted band size: 16 kDa
Observed band size: 16kDa