

MPO Polyclonal Antibody

(Catalog # A54284)

Description

MPO 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

P05164

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Myeloperoxidase protein (165-278AA)

Storage

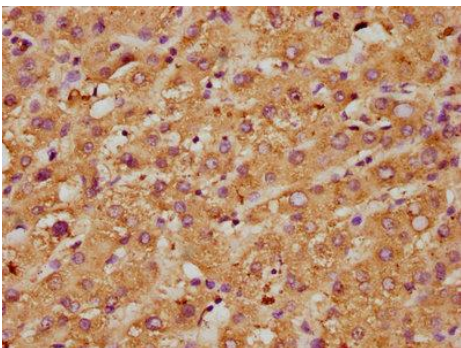
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

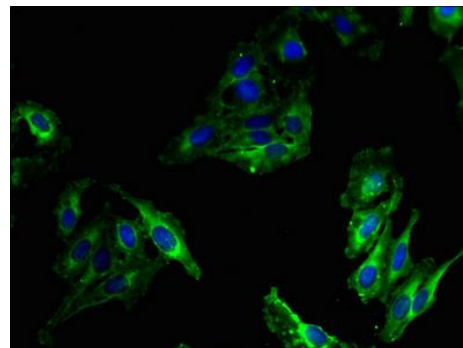
MPO

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:500-1:1000, IF:1:50-1:200



IHC image of MPO Antibody diluted at 1:600 and staining in paraffin-embedded human liver tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



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