

Nqo2 Polyclonal Antibody

(Catalog #A50973)

Background

The enzyme apparently serves as a quinone reductase in connection with conjugation reactions of hydroquinones involved in detoxification pathways as well as in biosynthetic processes such as the vitamin K-dependent gamma-carboxylation of glutamate residues in prothrombin synthesis.

Description

Nqo2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse, Rat

Isotype

IgG

Uniprot ID

Q6AY80

Purification

Protein G purified

Immunogen

Recombinant Rat Ribosyldihyronicotinamide dehydrogenase [quinone] protein (1-231AA)

Storage

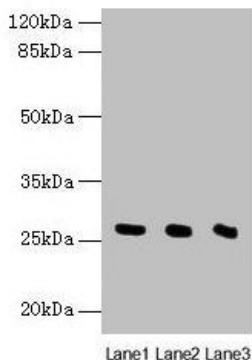
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

NRH dehydrogenase [quinone] 2, NRH:quinone oxidoreductase 2, Quinone reductase 2, QR2, Nqo2

Application

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



Western blot

All lanes: Nqo2 Polyclonal Antibody at 8µg/ml

Lane 1: Rat heart tissue

Lane 2: Mouse liver tissue

Lane 3: Mouse kidney tissue

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 27 kDa

Observed band size: 27 kDa