
DHRS3 Polyclonal Antibody, FITC Conjugated

(Catalog #A62412)

Background

Catalyzes the reduction of all-trans-retinal to all-trans-retinol in the presence of NADPH.

Description

DHRS3 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O75911

Purification

Affinity purified

Immunogen

Recombinant Human Short-chain dehydrogenase/reductase 3 protein (30-169AA)

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

Short-chain dehydrogenase/reductase 3, DD83.1, Retinal short-chain dehydrogenase/reductase 1, retSDR1, Retinol dehydrogenase 17, Short chain dehydrogenase/reductase family 16C member 1, DHRS3, RDH17, SDR16C1, UNQ2424/PRO4983

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
