

DAOA Polyclonal Antibody

(Catalog # A53792)

Description

DAOA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P59103

Purification

>95%, Protein G purified

Immunogen

Recombinant Human D-amino acid oxidase activator protein (1-153AA)

Storage

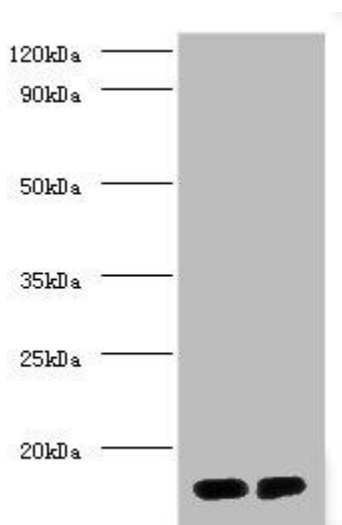
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein G72 DAOA G72

Application

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



Western blot
All lanes: DAOA Polyclonal Antibody at 2ug/ml
Lane 1: PC-3 whole cell lysate
Lane 2: mouse gonad tissue
Secondary
Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 19, 2, 10, 15 kDa
Observed band size: 18 kDa