

ANKRD23 Polyclonal Antibody

(Catalog # A65982)

Background

May be involved in the energy metabolism. Could be a molecular link between myofibrillar stretch-induced signaling pathways and muscle gene expression.

Description

ANKRD23 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

Q86SG2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Ankyrin repeat domain-containing protein 23 protein (1-200AA)

Storage

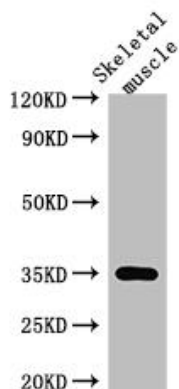
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

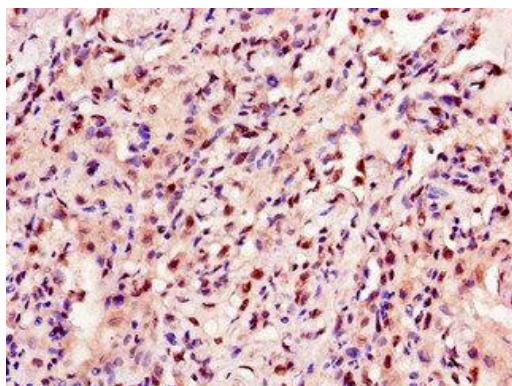
Ankyrin repeat domain-containing protein 23, Diabetes-related ankyrin repeat protein, Muscle ankyrin repeat protein 3, ANKRD23, DARP

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200



Positive WB detected in: Mouse skeletal muscle tissue
All lanes: ANKRD23 Polyclonal Antibody at 2.5 ug/ml
Secondary: Goat polyclonal to rabbit IgG at 1/50000 dilution
Predicted band size: 35,30 kDa; Observed band size: 35 kDa



Immunohistochemistry of paraffin-embedded human lung cancer using ANKRD23 Polyclonal Antibody at dilution of 1:100.