

LYPD2 Polyclonal Antibody

(Catalog # A65654)

Background

extracellular exosome

Description

LYPD2 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

Q6UXB3

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Ly6/PLAUR domain-containing protein 2 protein (23-103AA)

Storage

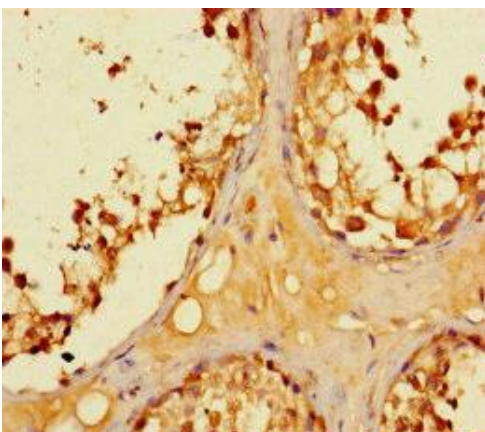
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

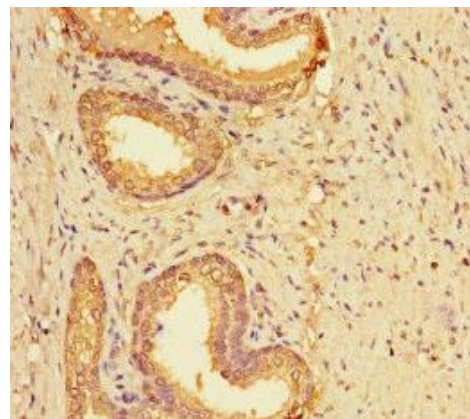
Ly6/PLAUR domain-containing protein 2, LYPD2, LYPDC2, UNQ430/PRO788

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human testis tissue using LYPD2 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human prostate cancer using LYPD2 Polyclonal Antibody at dilution of 1:100