

Mup11 Polyclonal Antibody

(Catalog # A54563)

Background

Binds pheromones that are released from drying urine of males. These pheromones affect the sexual behavior of females.

Description

Mup11 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

P04938

Purification

Protein G purified

Immunogen

Recombinant Mouse Major urinary protein 11 protein (1-151AA)

Storage

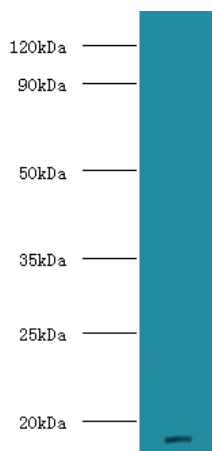
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

MUP11 and MUP8, Mup8, AND, Mup11

Application

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



Western blot

All lanes: Mup11 Polyclonal Antibody at 2ug/ml+mouse liver tissue

Secondary

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 18kDa

Observed band size: 18kDa