

RPS18 Polyclonal Antibody

(Catalog # A51757)

Background

Located at the top of the head of the 40S subunit, it contacts several helices of the 18S rRNA

Description

RPS18 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

P62269

Purification

>95%, Protein G purified

Immunogen

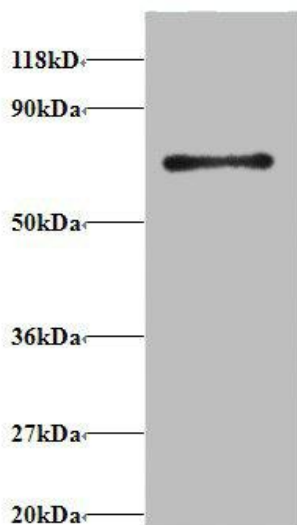
Recombinant Human 40S ribosomal protein S18 protein (2-152AA)

Storage

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

Application

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



Western blot
RPS18 Polyclonal Antibody at 2ug/ml+293T whole cell lysate
Secondary
Goat polyclonal to Rabbit IgG at 1/15000 dilution
Predicted band size: 16 kDa
Observed band size: 75 kDa