

SSB Polyclonal Antibody

(Catalog # A54756)

Background

Binds to the 3' poly(U) terminus of nascent RNA polymerase III transcripts, protecting them from exonuclease digestion and facilitating their folding and maturation.

Description

SSB 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

P05455

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Lupus La protein (1-408AA)

Storage

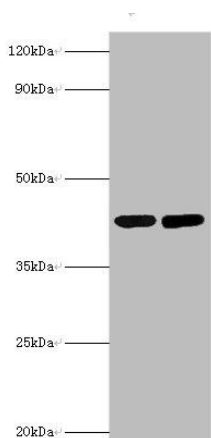
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

La autoantigen La ribonucleoprotein Sjogren syndrome type B antigen SSB

Application

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



Western blot
All lanes: SSB Polyclonal Antibody at 2ug/ml
Lane 1: mouse liver tissue
Lane 2: mouse brain tissue
Secondary
Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 47kDa
Observed band size: 47kDa