
SARS Polyclonal Antibody, HRP Conjugated

(Catalog # A53978)

Description

SARS Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P49591

Purification

Protein G purified

Immunogen

Recombinant Human Serine--tRNA ligase, cytoplasmic protein (2-233AA)

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

Serine--tRNA ligase, Short name=SerRS, Seryl-tRNA(Ser/Sec) synthetase, SARS, SERS

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
