





SARS glycoprotein M

SARS coronavirus glycoprotein M recombinant, *E. coli*

Cat. No.	Amount
PR-1453-S	200 µg
PR-1453-L	1 mg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 6 months

Molecular Weight: about 30.6 kDa

Form: liquid (Supplied in Tris-HCl 50 mM pH 8.0, Sodium Chlorid 60 mM, Glutathione 10 mM, Glycerol 50 %)

Applications:

Suitable for use in ELISA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not be excluded.

Specificity:

Immunoreactive with sera of SARS coronavirus-infected individuals

