



SARS-CoV-2 - nucleocapsid protein - 62.2 kDa

SARS-CoV-2 - coronavirus nucleocapsid protein
recombinant, *E. coli*

Cat. No.	Amount
PR-1456-S	200 µg
PR-1456-L	1 mg

For *in vitro* use only!

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Weight: about 62.2 kDa

Form: liquid (Supplied in Sodium Phosphate 100 mM, pH 8.0, Tris-HCl 210 mM, Sodium Chlorid 300 mM, Imidasol 500 mM, Urea 8 M)

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

This mixture contains dangerous substances



Signal word: Danger

Hazard statements

H315 Causes skin irritation.

H318 Causes serious eye damage.

H360D May damage the unborn child.

Precautionary statements

P201 Obtain special instructions before use.

P302 + P352 IF ON SKIN: Wash with plenty of water.

P305 + P351 + P338 IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

P308 + P313 IF exposed or concerned: Get medical advice/attention.

P332 + P313 If skin irritation occurs: Get medical advice/attention.

P362 + P364 Take off contaminated clothing and wash it before reuse.

For further information see Material Safety Data Sheet.