

Covid-19 Nucleocapsid Antibody

Cat. No. C15310269

Type: Polyclonal	Specificity: Covid-19
Size: 100 µl	Isotype: NA
Concentration: not determined	Host: Rabbit
Lot No.: A3259-002	Purity: Whole antiserum from rabbit containing 0.05% azide.
Storage buffer: NA	Storage conditions: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Last Data Sheet Update: August 12, 2020

Description

Polyclonal antibody raised in rabbit against the structural nucleocapsid (NC) protein from the Covid-19 virus using a recombinant protein.

Applications

Applications	Suggested dilution	References
Western Blotting	1:1,000	Fig 1

Target Description

The nucleocapsid protein is one of the 4 structural proteins of the Covid-19 virus. This protein has non-specific nucleic acid binding capacities and associates with the viral RNA. It facilitates viral replication and is involved in viral particle assembly and release from the cell. The NC protein is a highly immunogenic and abundantly expressed protein during infection and is a frequently used target in vaccine development and serological assays.

Validation data

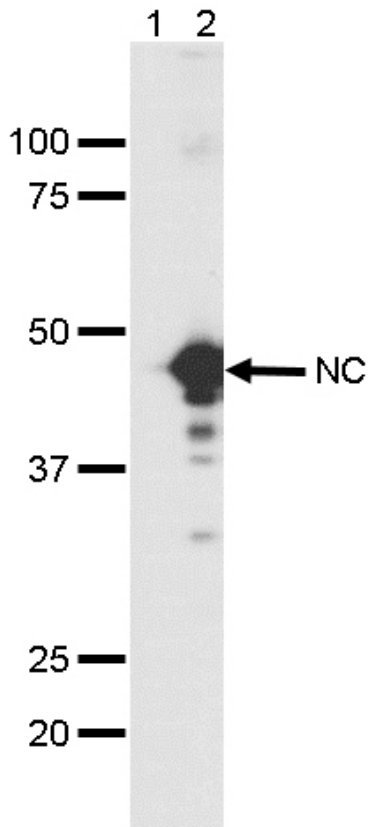


Figure 1. Western blot analysis using the Diagenode antibody directed against Covid-19 NC

Western blot was performed on 2 μ g of purified Covid-19 nucleocapsid (NC) protein (6xHis tagged) using the Diagenode antibody against Covid-19 NC (Cat. No. C15310269), diluted 1:1,000 in PBS-T containing 3% NFDM. The marker is shown on the left, the position of the NC protein is indicated on the right. Lane 1 shows the Western blot analysis with the pre-immune serum, used as a negative control.