

## Covid-19 Spike Antibody

**Cat. No. C15310268**

Type: Polyclonal	Specificity: Covid-19
Size: 100 µl	Isotype: NA
Concentration: not determined	Host: Rabbit
Lot No.: A3257-002	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 Spike protein from the Covid-19 virus using a recombinant protein.

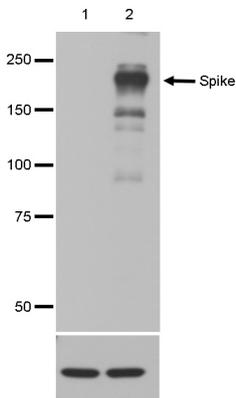
### Applications

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

### Target Description

Spike is one of the 4 structural proteins of the Covid-19 virus. It's a trimeric transmembrane glycoprotein precursor that contains 1273 amino.acids and is cleaved into two segments S1 and S2. The spike protein covers the surface of the virus and is responsible for binding to the ACE2 receptor and entry into the host cell. It is therefore to most important target for the generation of a vaccine.

**Validation data**



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

Western blot was performed on protein extracts from HEK293 cells transfected with Spike-P2A-GFP (lane 2) and on untransfected cells (lane 1) using the Diagenode antibody against Covid-19 Spike (cat. No. C15310268), diluted 1:1,000 in PBS-T containing 3% NFD. The marker is shown on the left, the position of the Spike protein is indicated on the right.