

AbFlex® SARS-CoV-2 Spike Antibody (rAb) (AM002414)

Catalog Nos: 91363, 91364

RRID: AB_3216349

Clone: AM002414 (414-2)

Application(s): ELISA, NEU

Reactivity: Virus

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Isotype: IgG2a

Molecular Weight: 141 kDa (full length S1 protein)

Background: AbFlex® antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports. AbFlex® SARS-CoV-2 Spike (rAb) (AM002414) antibody was expressed in was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

COVID-19, which is short for coronavirus disease 2019, is the official name of the respiratory disease caused by infection with the novel coronavirus SARS-CoV-2. The virus that causes COVID-19 was named SARS-CoV-2 because it is a coronavirus genetically similar to, yet distinct from, the virus that caused the severe acute respiratory syndrome (SARS) outbreak in 2003. Studying the details of how this virus replicates and causes the disease will allow scientists and physicians to more rapidly develop fast and accurate methods of detection as well as to deploy therapeutic and vaccine strategies.

Immunogen: N/A - derived from COVID-19 patients who have cleared the virus. Their antibodies were screened for reactivity to SARS-CoV-2 and then sequenced and expressed as recombinant antibodies.

Buffer: 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, and 30% glycerol. Sodium azide is highly toxic.

Application Notes:

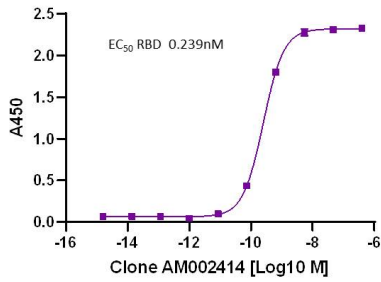
Applications Validated by Active Motif:

ELISA: 0.05 – 1 µg/ml

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

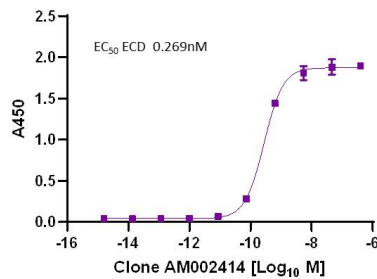
SARS-CoV2 Spike AbFlex Antibody
Clone AM002414 EC₅₀ with RBD



AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM002414) tested by ELISA using SARS Spike protein RBD.

SARS-CoV-2 Spike RBD protein was coated onto microtiter plates at 10 µg/mL and then incubated with a dilution series of AbFlex SARS-CoV-2 Spike Antibody (rAb) (clone AM002414). Bound antibodies were detected with anti-mouse IgG conjugated to horseradish peroxidase (HRP) followed by incubation with HRP Substrate and then measuring the resulting absorbance at 450 nm.

SARS-CoV2 Spike AbFlex Antibody
Clone AM002414 EC₅₀ with ECD



AbFlex SARS-CoV-2 Spike Antibody (rAb) (AM002414) tested by ELISA using SARS Spike protein ECD.

SARS-CoV-2 Spike Extracellular Domain (ECD a.a. 14-1213) protein was coated onto microtiter plates at 5 µg/mL and then incubated with a dilution series of AbFlex SARS-CoV-2 Spike Antibody (rAb) (clone AM002414). Bound antibodies were detected with anti-mouse IgG conjugated to horseradish peroxidase (HRP) followed by incubation with HRP Substrate and then measuring the resulting absorbance at 450 nm.