

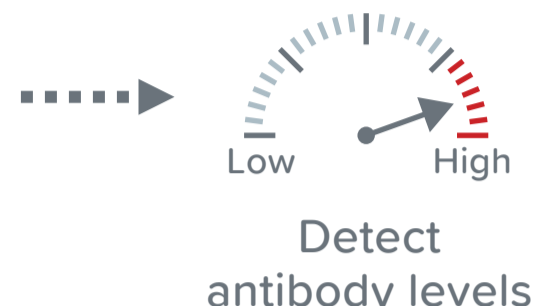
DETECTING SARS-COV-2 ANTIBODIES IN SERUM AND PLASMA SAMPLES

Convalescent plasma therapy is a potential treatment under investigation where antibodies from recovered patients are transfused to current COVID-19 patients with the intent to help them fight the infection and buy time until their immune system can produce antibodies. However, not all recovered persons have the same amount of antibody titers suitable for such transfusions. Reliable, simple, and rapid methods are needed to assess relative antibody titers in plasma samples.

HOW WILL YOU KNOW IF PLASMA ANTIBODY LEVELS ARE SUFFICIENT?

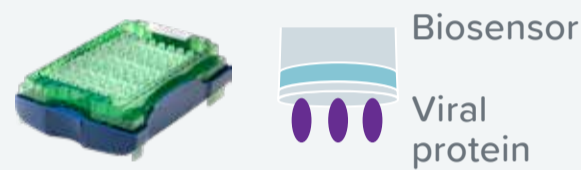


.....
Plasma separation from blood



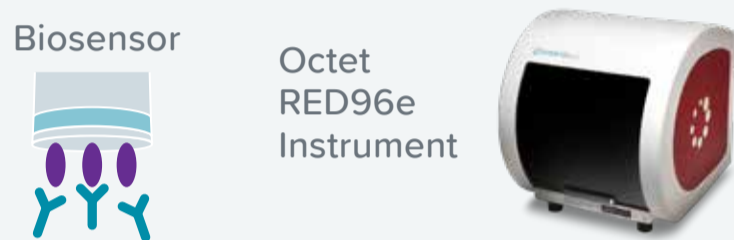
HOW TO DETECT PLASMA ANTIBODIES USING THE OCTET SYSTEM

Step 1: Prepare a coronavirus antibody specific biosensor



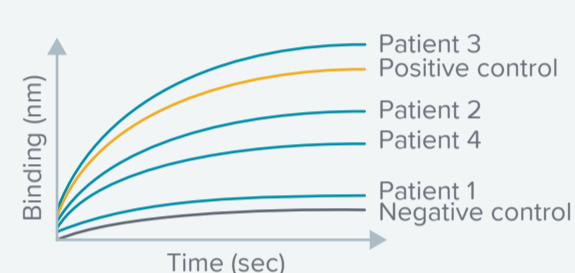
Immobilize a virus protein such as the receptor binding domain (RBD) of the SARS CoV-2 spike protein.

Step 2: Detect antibody binding signals



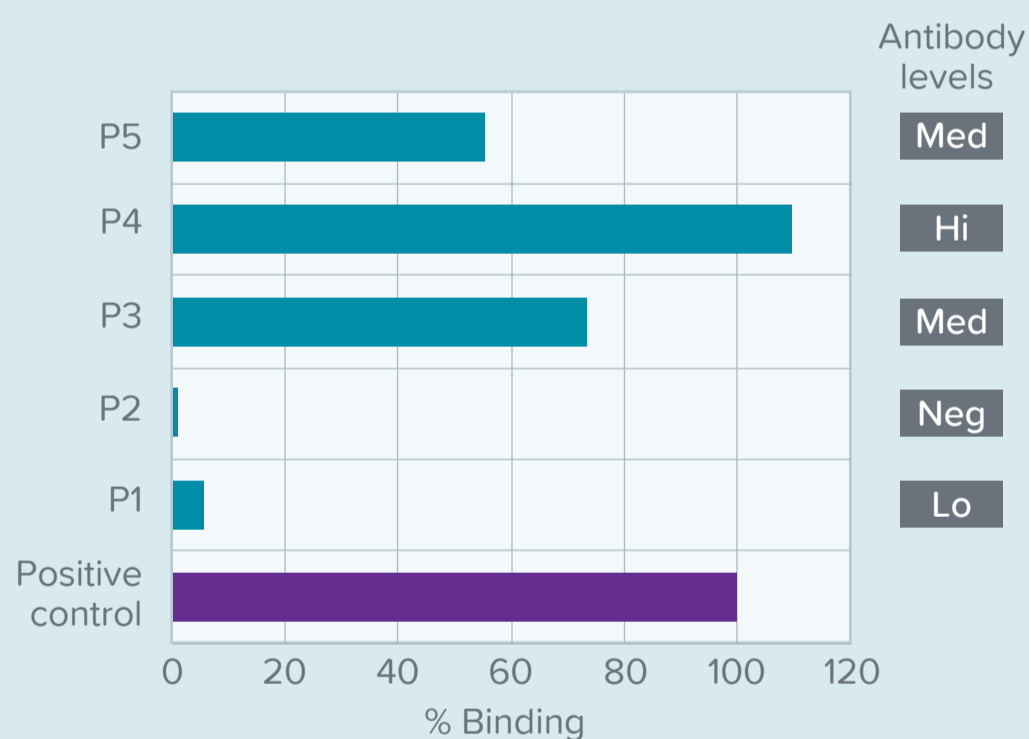
Dip the coronavirus biosensor into diluted patient plasma or serum samples. Block the biosensor with non-relevant serum or blocking buffer if needed to prevent non-specific binding.

Step 3: Analyze binding data



Quantitate relative antibody levels in plasma or serum.

WHAT THE DATA TELLS YOU



HOW DO OCTET SYSTEMS HELP YOU GET DATA FAST?

- Rapid test generating real-time data
- Simple workflow and easy to use
- Low cost assay development
- Wide-range of available biosensor chemistries to immobilize virus proteins

A RANGE OF BLI PLATFORMS TO MEET YOUR NEEDS

Octet HTX Instrument

- 96-channel fully automated
- Samples in 96/384-well plates
- Robot compatible



Octet RED384 Instrument

- 16-channel fully automated
- Samples in 96/384-well plates
- Robot compatible



Octet RED96e Instrument

- 8-channel fully automated
- Samples in 96-well plates



Octet K2 Instrument

- 2-channel fully automated
- Samples in 96-well plates



BLitz Instrument

- Single channel manual system
- Sample consumption from 4–250 µL



Learn more about how Octet systems are used in COVID-19 research at www.fortebio.com/covid19research