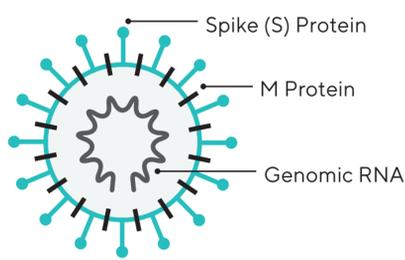


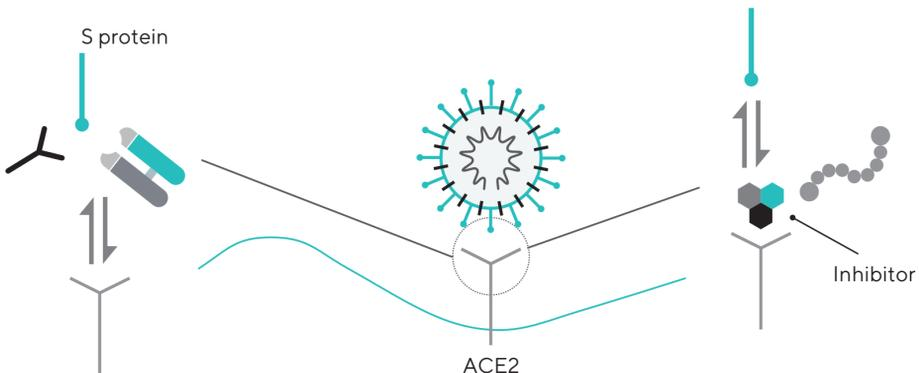
# Understanding COVID-19 Vaccine and Therapeutic Development Research

## SARS-CoV-2 Mechanism for Infecting Hosts

Coronavirus uses its spike (S) protein to attach onto ACE2 receptors on human cells. Targeting this virus-host receptor interaction can prevent infection.



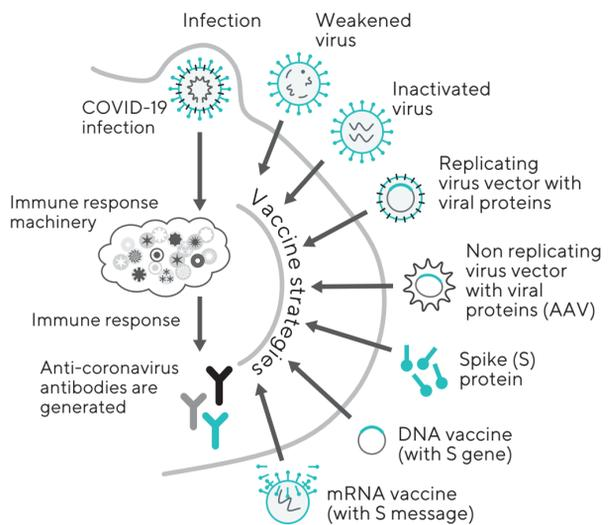
## Targeting Virus-Host Receptor Recognition to Block Viral Entry



- Developing antibodies, FAB, and scFv molecules that bind to viral spike proteins are at the forefront at inhibiting viral entry. Similarly, small-molecule and peptides that target the spike protein are also evaluated as alternate therapeutic strategies.
- The selection and characterization of lead candidates based on accurate target binding kinetics, affinities and inhibitory potency is vital in the discovery workflow.

## Strategies for COVID-19 Vaccine Development

The scientific community has made tremendous efforts to understand the disease, and unparalleled efforts are ongoing to develop vaccines and treatments. Studying antibodies generated from COVID-19 infections and vaccines can provide invaluable information towards therapeutic antibody development and engineering. Kinetics, affinity and epitope interrelationships are important factors in vaccine and therapeutic design.

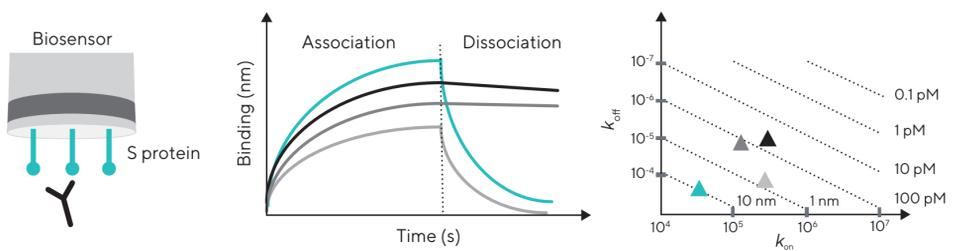


## Octet® Systems: All-In-One Platform for Therapeutic Investigation and Bioproduction

Octet® systems perform critical measurements with the speed and reliability needed for investigating therapeutic strategies and producing novel medicine.

## Evaluate Binding Kinetics and Affinities in Real-Time

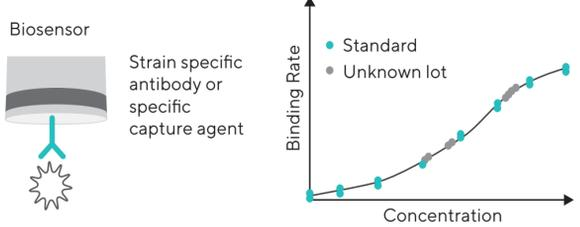
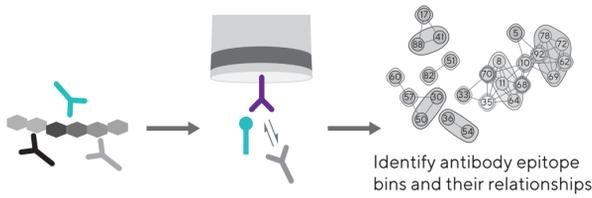
Pick lead candidates with the best kinetics and affinity properties



## Assess Epitope Diversity and Coverage for Greater Success

## Perform Titer, Potency and Stability Measurements

Epitope binning assays help identify antibodies that block the same epitope on a target antigen and are crucial when it comes to identifying or engineering mAbs with favorable kinetics and affinity profiles.



- Three epitope binning assay formats to choose between.
- Different Bio-Layer Interferometry (BLI) systems to meet your throughput needs and budget.

- Fast at-line testing – Quantitate a 96-well plate for AAV titers within minutes.
- Robust assays for antibody and virus titer (ex. Influenza, AAV), potency and stability determination in upstream and downstream bioprocess development.

## How Do Octet® Systems Help You Get Data Fast?

- Simple Dip and Read workflows help you develop assays quickly and get results within minutes.
- Minimal sample preparation time with analysis possible directly from crude samples.
- Range of available biosensor chemistries makes assay development easy.

Learn more about how Octet® systems are used in COVID-19 research at [sartorius.com/covid19](https://www.sartorius.com/covid19)