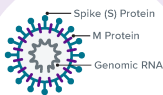


# UNDERSTANDING COVID-19 VACCINE AND THERAPEUTIC DEVELOPMENT RESEARCH

## 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.



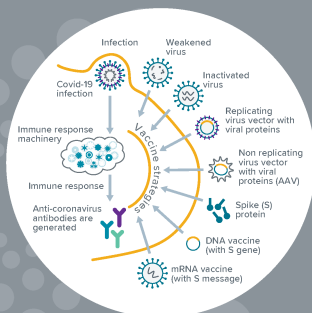
## TARGETING VIRUS-HOST RECEPTOR RECOGNITION TO BLOCK VIRAL ENTRY



- 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 affinities, kinetics and inhibitory potency is vital in the discovery workflow.

## CURRENT STRATEGIES FOR COVID-19 VACCINE DEVELOPMENT

Studying antibodies generated from Covid-19 infections and vaccines can provide invaluable information towards therapeutic antibody development and engineering. Affinity, kinetics, and epitope interrelationships are important factors in vaccine and therapeutic design.

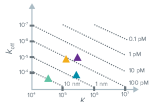
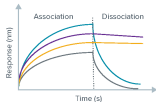


## OCTET: ALL-IN-ONE PLATFORM FOR THERAPEUTIC INVESTIGATION AND BIOPRODUCTION

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

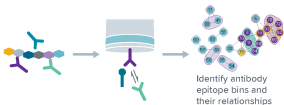
## EVALUATE BINDING KINETICS AND AFFINITIES IN REAL TIME

Pick lead candidates with the best kinetic and affinity properties



## ASSESS EPIOTOPE DIVERSITY AND COVERAGE FOR GREATER SUCCESS

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.

## PERFORM TITER, POTENCY AND STABILITY MEASUREMENTS

Biosensor



- 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 construction easy.

Learn more about how Octet systems are used in COVID-19 research at [www.fortebio.com/covid19research](http://www.fortebio.com/covid19research)

**FORTÉBIO**  
A SARTORIUS BRAND