

Data Sheet

Antibody Neutralization Assay with SARS-CoV-2-614G Pseudovirus Particles (SARS-CoV-2-614G-PP)

Features

- **Robust:** Excellent signal to noise (basal) ratio.
- **Easy to use:** Amenable to HTS format (96-well, 384-well and 1536-well format).

Applications

- Working perfectly for Luc Pseudovirus to get robust signal, screening potential neutralization antibodies against SARS-CoV-2

Product Information

Catalog Number:

Components

CB-97100-146G-1ml**1ml SARS-CoV-2-614G Pseudovirus Particles****CB-97100-146G-5ml****5ml SARS-CoV-2-614G Pseudovirus Particles****Storage**

Store at -80°C

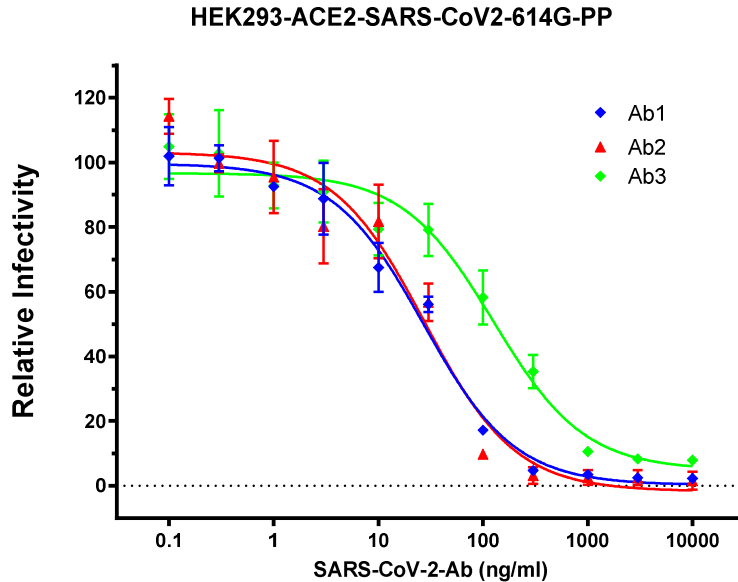
ASSAY PROTOCOL**Cell Infection:**

1. Count HEK293-ACE2 cells (CB-97100-203) to be infected and seed ~7.5K cells per well into 384-well plates (15 µl per well) in DMEM with 10% HyClone™ FetalClone™ II Serum (no antibiotics).
2. On the 2nd day, dilute the antibodies (5X of the final concentrations) in DMEM with 10% HyClone™ FetalClone™ II Serum (no antibiotics).
3. Take 5 µl of the antibody solution. Add into the well containing 12.5 µl of SARS-CoV-2-PP.
4. Incubate at 37°C for 1 hr.
5. remove media, add 17.5 µl of SARS-CoV-2-PP and antibody mix into each well. Centrifuge at 700 rpm for 15 min at 4°C
6. Add additional 7.5 µl of DMEM with 10% FC into each well.
7. Incubate at 37 °C for 42 hr

Measurement of Luciferase Activity in Infected cells

1. Remove supernatant
2. Add 20 µl Codex's Luciferase assay reagent (CB-80552-010) into each well.
3. Read in a luminescence plate reader, record the data.

DATA



	Ab1	Ab2	Ab3
Bottom	0.2185	-1.720	4.886
Top	99.66	103.2	96.69
LogIC50	1.432	1.439	2.099
IC50	27.02	27.50	125.7
Span	99.45	104.9	91.80

Figure 1. Antibody Neutralization Assays with SARS-CoV-2-614G PP