

## Data Sheet

## Codex SARS-CoV-2 Pseudovirus Particles (SARS-CoV-2-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 inhibitor to block SARS-CoV entry and viral protein translation

**Product Information**

Catalog Number:

Components

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

Store at -80°C

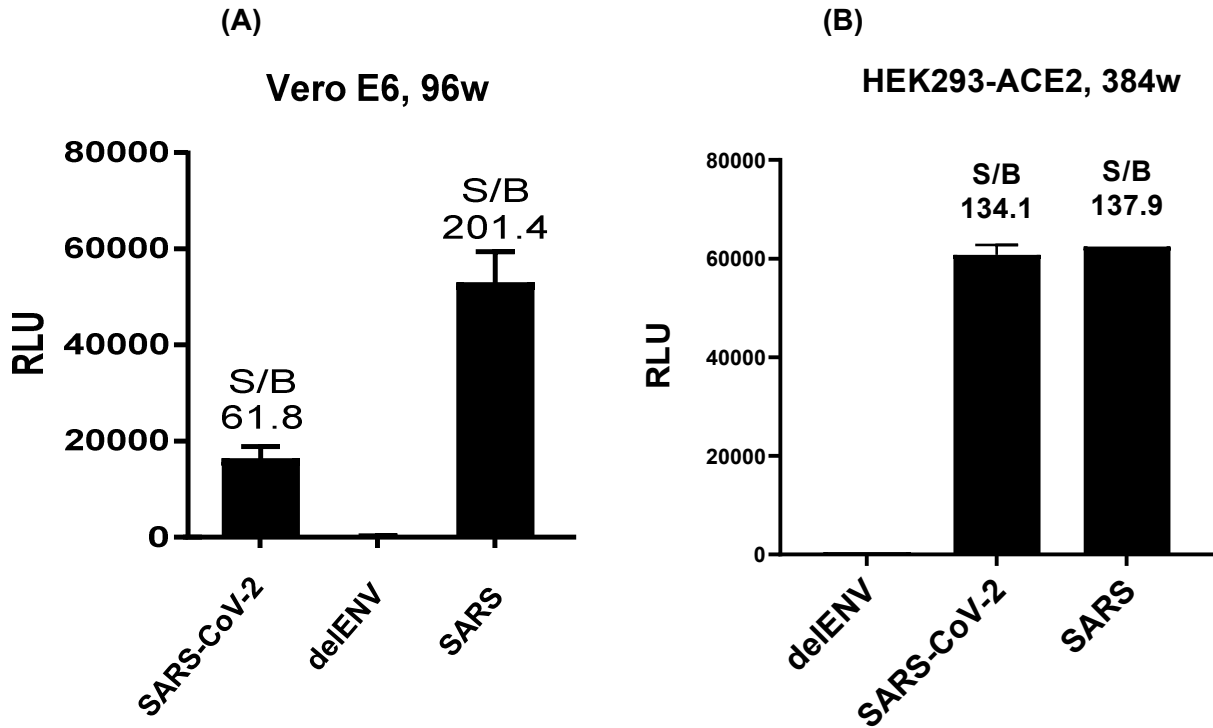
**ASSAY PROTOCOL****Cell Infection:**

1. Count Vero E6 cells/HEK293-ACE2-GFP cells (CB-97100-203) to be infected and seed ~20K cells per well into 96-well plates (50 µl per well) DMEM with 10% FC (no antibiotics) or 5K cells per well into 384-well plates (15 µl per well)
2. Culture cells overnight to make sure the cells stably adhere to the plates.
3. On the 2<sup>nd</sup> day, remove media, add 50 µl SARS-CoV-2-PP into each well (12.5 µl for 384-well plate). Spin at 700 rpm for 15 min at 4°C
4. Incubate for 2 hr at 37 °C
5. Add 50 µl DMEM with 10% FC into each well (12.5 µl for 384-well plates).
6. Incubate for 48 hr at 37 °C

**Measurement of Luciferase Activity in Infected cells**

1. Remove supernatant
2. Add 100 µl Codex's Luciferase assay reagent (CB-80552-010). (20 µl for 384-well plates).
3. Read in a luminescence plate reader, record the data.

DATA



**Figure 1. Pseudoviral Particle (PP) Infection Assays**

(A) SARS and SARS-CoV-2 pseudoviral particles on Vero E6 cells in 96-well format.

(B) SARS and SARS-CoV-2 pseudoviral particles on HEK293-ACE2 cells in 384-well format.

**Legends:** **SARS-CoV-2:** SARS-CoV-2 MLV Pseudovirus Particles;  
**SARS:** SARS MLV Pseudovirus Particles.  
**deIENV:** MLV Pseudovirus Particles without Spike protein