

## Data Sheet

# Codex SARS-CoV-2-XBB.1.16.6 Variant Pseudovirus Particles (SARS-CoV-2-XBB.1.16.6-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-2 XBB.1.16.6 variant entry and viral protein translation

**Product Information**

Catalog Number:

Components

CB-97200-213-1ml

1ml SARS-CoV-2-XBB.1.16.6 Variant Pseudovirus Particles

CB-97200-213-5ml

5ml SARS-CoV-2- XBB.1.16.6 Variant Pseudovirus Particles

**Storage**

Store at -80°C

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

1. HEK293-ACE2 cells (CB-97100-203) to be infected and seed ~20K cells per well into 96-well plates (50 µl per well) DMEM with 10% HyClone™ FetalClone™ II Serum (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-XBB.1.16.6-PP into each well (12.5 µl for 384-well plate). Spin at 700 rpm for 15 min at 4°C

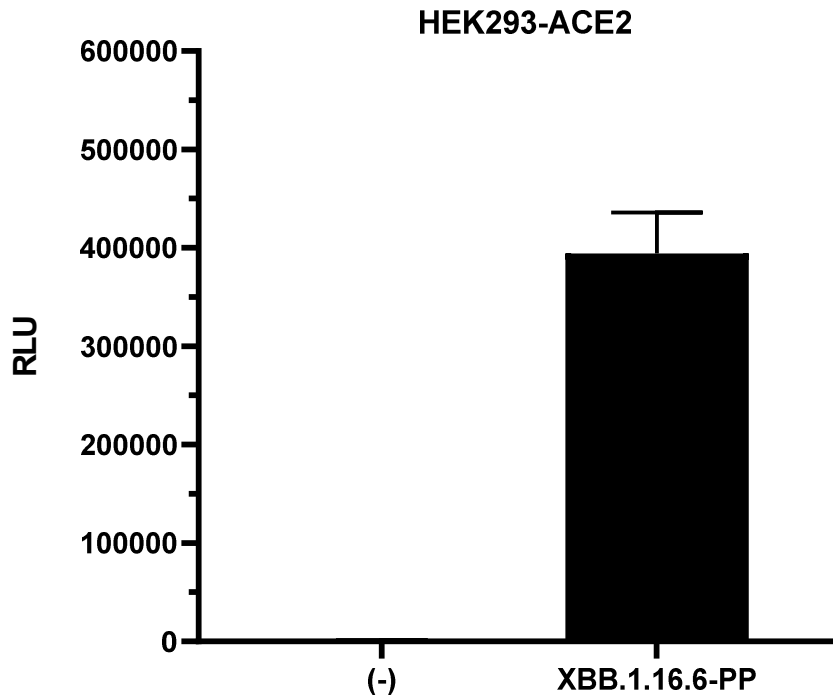
**Note: Thaw the PPs quickly with room temperature water**

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 42 hr at 37 °C

**Measurement of Luciferase Activity in Infected cells**

1. Add 100 µl Codex's Luciferase assay reagent (CB-80552-010). (25 µl for 384-well plates).
2. Read in a luminescence plate reader, record the data.

## DATA



**Figure 1. Pseudoviral Particle (PP) Infection Assays.** SARS-CoV-2-XBB.1.16.6 variant pseudoviral particles on HEK293-ACE2 cells in 384-well format.

**Legends:** SARS-CoV-2-XBB.1.16.6-PP: SARS-CoV-2-XBB.1.16.6 Variant MLV Pseudovirus Particles;

**Figure 2. Antibody neutralization assays using XBB.1.16.6 pseudoviral particle (PP) infection method.**

Each antibody dose-response curve represents the relative infectivity under serial dilutions of antibody. At each Ab concentration, HEK293-ACE2 cells were co-incubated with SARS-CoV-2 XBB.1.16.6 pseudoviral particles in the presence of the antibody.