

Data Sheet

Codex SARS-CoV-2-DELTA Variant Pseudovirus Particles (SARS-CoV-2-DELTA-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 DELTA variant (B.1.167.2) entry and viral protein translation

Product Information

Catalog Number:

Components

CB-97100-161-1ml**1ml SARS-CoV-2-DELTA Variant Pseudovirus Particles****CB-97100-161-5ml****5ml SARS-CoV-2-DELTA 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 2nd day, remove media, add 50 µl SARS-CoV-2-DELTA-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 µ 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. 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

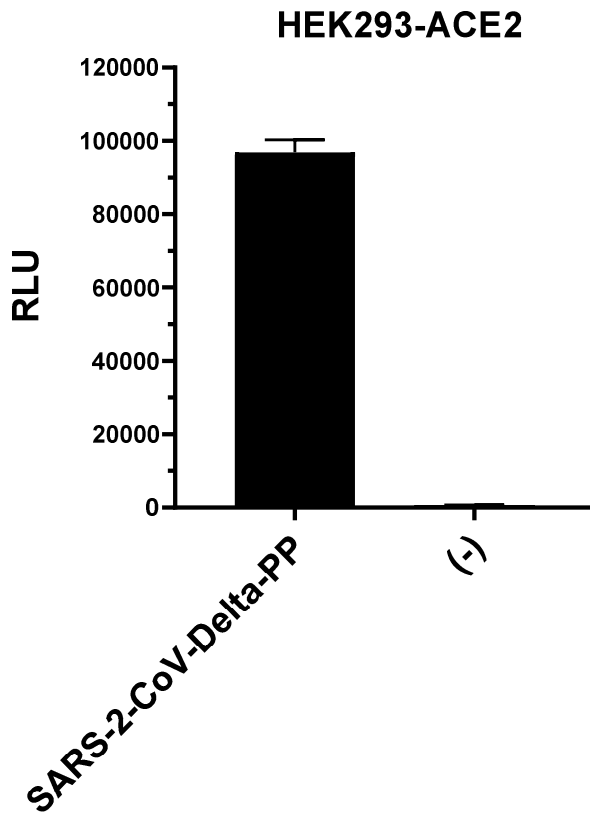


Figure 1. Pseudoviral Particle (PP) Infection Assays

SARS-CoV-2-DELTA variant pseudoviral particles on HEK293-ACE2 cells in 96-well format.

Legends: SARS-CoV-2-DELTA-PP: SARS-CoV-2-DELTA Variant MLV Pseudovirus Particles;