

# Genedrive<sup>®</sup> 96 SARS-CoV-2 Kit Quick Reference Guide



## Step 1

Place 96 well plate in a cool block **for the duration of reaction setup.**



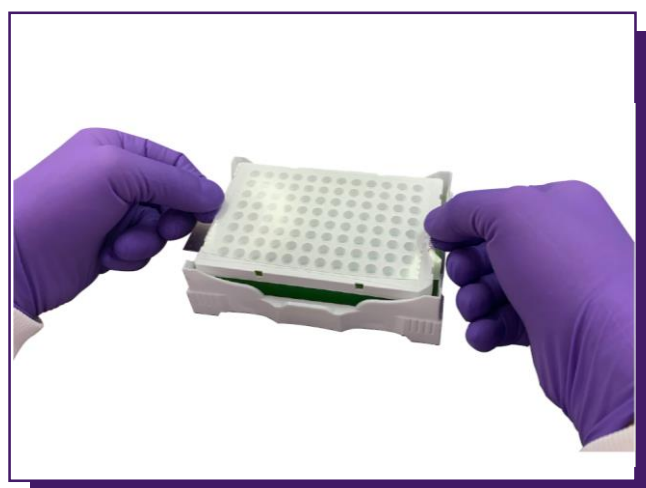
## Step 2

Carefully peel back the foil seal. It is recommended to do this in sections to minimise open wells during pipetting. Check there is a reagent bead in each well.



## Step 3

Pipette 20µL of extracted specimen or control into the side of each well. Add **20µL of nuclease free water to any wells that do not contain a specimen. Ensure all beads have been rehydrated**



## Step 4

Carefully peel optical film seal from its backing and seal the plate. Use of a microtitre sealing brayer or similar is recommended.



## Step 5

Carefully drop cool block 3 times from 1 cm height. Vortex the cool block/plate for **3 x 15 seconds at full power (total 45 seconds).**



## Step 6

Carefully drop cool block 3 times from 1 cm height. Incubate on cool block for **5 minutes.**



## Step 7

Vortex the cool block/plate for **3 x 15 seconds at full power (total 45 seconds).**



## Step 8

Centrifuge the plate **2 minutes at 2000 RPM.**



## Step 9

Load the plate on the Real-Time PCR instrument and start the program immediately. Ensure all the run parameters are correct before starting.