

Genedrive® 96 SARS-CoV-2 Kit

The Genedrive® 96 SARS-CoV-2 Kit is a **unique, one-step test** utilising a single PCR bead format provided in a 96 well plate **adaptable for automation**, which makes it an ideal solution to **increase throughput in your laboratory**.

In external evaluations the Genedrive® 96 SARS-CoV-2 Kit achieved a sensitivity of 100% and a specificity of 99%. The dual gene target design gives greater inclusivity than tests with a single gene target.

Validated platforms

- Validated on multiple market leading Real-Time PCR platforms for high throughput testing to meet the needs of larger laboratories and high volume in the COVID-19 pandemic
 - ABI 7500 Fast
 - Roche LightCycler® 480 II
 - Bio-Rad CFX96

Fast result analysis

- Free Exporter software for instant result analysis

Relevant gene targets

- Unaffected by recently reported new Covid-19 variants* from UK, South Africa and Brazil

Ease of use

- Single PCR bead format – eliminates need for reagent preparation
 - Load extracted specimen and run the test
 - Up to a 50% reduction in the use of tips and 96 well plates than other RT-PCR kits
 - Automation compatible 96 well design

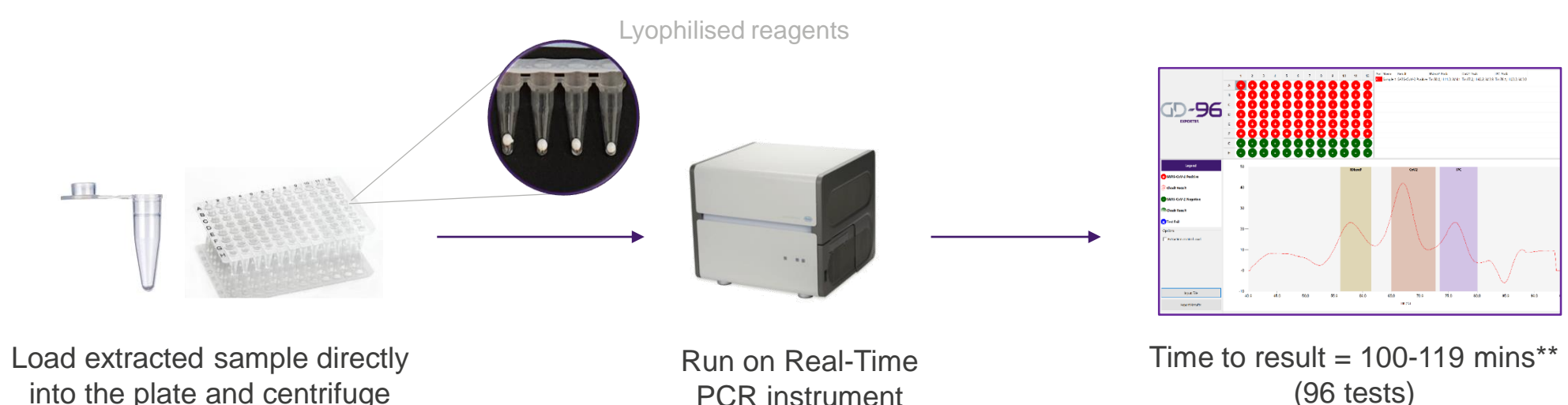
Excellent performance

- Sensitivity = 100%
- Specificity = 99%

Ambient storage

- Lyophilised reagents for ease of storage

Simple one-step procedure



For further information contact us at: enquiries@genedrive.com