

Covid-19 GeoMx-formatted Panel Probe R Mix ab273595

Overview

Product name Covid-19 GeoMx-formatted Panel Probe R Mix

Product overview Covid-19 GeoMx-formatted Panel Probe R Mix ab273595 is used with Covid-19 GeoMx-formatted Antibody Panel **ab273594**.

It contains:

- ID0314 - Probe R_ 9.2
- ID0315 - Probe R_ 9.3
- ID0321 - Probe R_ 9.4
- ID0322 - Probe R_ 9.5
- ID0318 - Probe R_ 9.6

IMPORTANT TECHNICAL NOTES:

Storage instructions: Shipped on dry ice. Store at -20°C long term.

Storage buffer: 120 µl in TE Buffer pH 8.0 (10mM TrisHCl, 0.1mM EDTA)

Concentration: 5nM for each Probe

Use this Probe R as explained in the GeoMx DSP nCounter Readout User Manual from NanoString (www.nanostring.com/geomx-online-user-manual). Each 120 µl of Probe R Mix is sufficient for 12 slides at 48 regions of interest per slide, or 576 total regions of interest.

For a full list of pre-validated NanoString GeoMx Assays, visit <http://www.nanostring.com/products/geomx-digital-spatial-profiler/geomx-assays>

To learn more about Abcam's GeoMx-formatted antibodies and custom conjugation services, please visit **here**.

NanoString, nCounter and GeoMx are trademarks or registered trademarks of NanoString Technologies, Inc., in the United States and/or other countries.

Properties

Storage instructions

Store at -80°C. Please refer to protocols.

Components	120 µl
Probe R Mix	1 x 120µl

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors