

Important Note

Significant updates to *therascreen*® PIK3CA RGQ PCR Kit Instructions for Use

Dear valued *therascreen* customer,

This is to inform you that we recently made some significant changes to the *therascreen* PIK3CA RGQ PCR Kit Instructions for Use (HB-2614). Updates to the IFU (R2) detail the steps taken to demonstrate the repeatability and validity of Q546R mutation calling and required updates to Q546R mutation calling/results demonstration that support the inclusion of identified Q546R mutations.

Key edits to the IFU

- **Updates to *therascreen* PIK3CA RGQ PCR Kit assay profiles**

The *therascreen* PIK3CA FFPE and Plasma Assay Profiles have been updated to address the increased probability of false Q546R mutation positive results caused by non-specific molecular interactions within the Q546R reaction and reinstate accurate detection of Q546R mutations positive results. Consequently, the Important Note for *therascreen* PIK3CA RGQ PCR Kit (HB-2854-001) dated December 2020 **no longer applies**.

Note: The updated *therascreen* PIK3CA FFPE and Plasma assay profiles are available for download from the related product page at www.qiagen.com. Go to **Resources** > **Protocols** to download the applicable assay profile.

Customers must update to these latest versions of both FFPE and Plasma assay profiles before reinstating use of Q546R calls:

- **therascreen_PIK3CA_FFPE_MDx Assay Profile Version 1.0.1:** For analysis of tissue specimens.
- **therascreen_PIK3CA_Plasma_MDx Assay Profile Version 1.0.2:** For analysis of plasma specimens.

The *therascreen* PIK3CA RGQ PCR Kit is specifically designed for use with Rotor-Gene® Q MDx (US) instrument operating with a computer installed with:

- Rotor-Gene Q AssayManager® v2.1
- Gamma MDx Plug-in version 1.0.0
- *therascreen_PIK3CA_FFPE_MDx* Assay Profile version 1.0.1 for analysis of tissue specimens
- *therascreen_PIK3CA_Plasma_MDx* Assay Profile version 1.0.2 for analysis of plasma specimens

Refer to the *Rotor-Gene Q MDx (US) User Manual* information regarding the Rotor-Gene Q MDx (US) instrument. The instrument must be maintained according to the requirements in the user manual.

Refer to *Rotor-Gene AssayManager v2.1 MDx Core Application User Manual* and *Rotor-Gene AssayManager v2.1 Gamma MDx (US) Plug-in User Manual* for further information regarding the software.

- **Additional clinical and performance data**

- Updated repeatability and reproducibility data for samples containing mutations in Q546R, including C420R, E542K, E545A, E545D, E545G, E545K, Q546E, H1047L, H1047R, and H1047Y
- Revised Clinical performance section to update progression-free survival (PFS) analysis, concordance, and participant/sample testing data following reanalysis with updated software parameters

- **Troubleshooting and Interpretation**

- Enhanced troubleshooting guidance for invalid sample flags
- Added clarification on cross-reactivity handling in MAF samples

- **Specimen Handling and Stability**

- Clarified specimen storage conditions and maximum storage times for FFPE tissue and plasma
- Update to General precautions specifically relating to guidance on handling and disposal of *Taq* DNA polymerase

For full details, refer to the latest version of the Instructions for Use available at www.qiagen.com

If you have any questions or concerns, please contact your local QIAGEN Technical Service Department through any of the following:

Telephone: 800 362 7737

Email: TechService-NA@qiagen.com

Best regards,
QIAGEN