

# Performance Evaluation of a Highly Mutation-Resilient RT-PCR Assay to Address Diagnostic Challenges with SARS-CoV-2 Variants of Concern

Jasmin Köffer<sup>1</sup>, Zoe Pounce<sup>2</sup>, Tammy Pham<sup>2</sup>, Emily Hodgson<sup>2</sup>, Xuemei Wang<sup>2</sup>, Géraldine Mercier<sup>2</sup>, Jelena Feenstra<sup>2</sup>, Manoj Gandhi<sup>2</sup>, Ulrich Eigner<sup>1</sup>

<sup>1</sup>MVZ Labor Dr. Limbach & Kollegen GbR, Heidelberg, Germany; <sup>2</sup>Thermo Fisher Scientific, South San Francisco, CA, USA

## Introduction

One of the considerations regarding SARS-CoV-2 variants of concern (VOC) is the impact of mutations on the performance of diagnostic tests. The TaqPath™ COVID-19 RNase P Combo Kit 2.0\* was designed with a total of 8 targets in 3 genes: *orf1a*, *orf1b* and *N* to compensate for emerging SARS-CoV-2 variants. In the current study we evaluated the clinical performance of the TaqPath™ COVID-19 RNase P Combo Kit 2.0\* as compared to the cobas® SARS-CoV-2 Assay (2 targets, 1 each in *orf1ab* and *E*-gene).

## Methods

The retrospective study was performed on 675 nasopharyngeal swab samples collected in Germany from February to August 2021. SARS-CoV-2 status was characterized using the cobas® SARS-CoV-2 Assay. The samples were tested using both tests in a blinded, randomized fashion and Positive Percent Agreement (PPA) and Negative Percent Agreement (NPA) were assessed. Discordant results were resolved using the TaqPath™ COVID-19 RT-PCR CE-IVD kit\* and sensitivity and specificity were calculated. Variant status for 116 samples was determined using the TIB MOLBIOL SARS-CoV-2 combi assay (Figure 1).

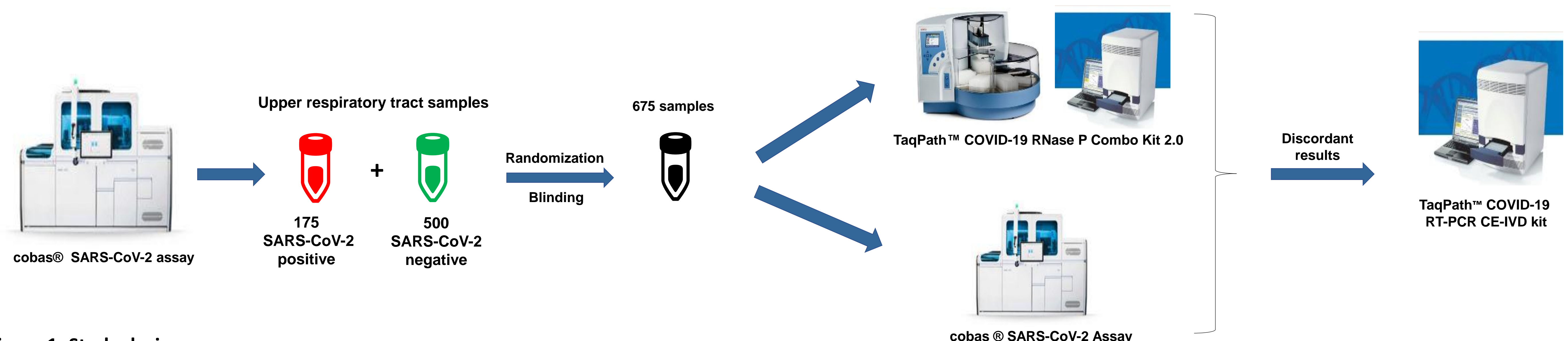


Figure 1: Study design

## Results

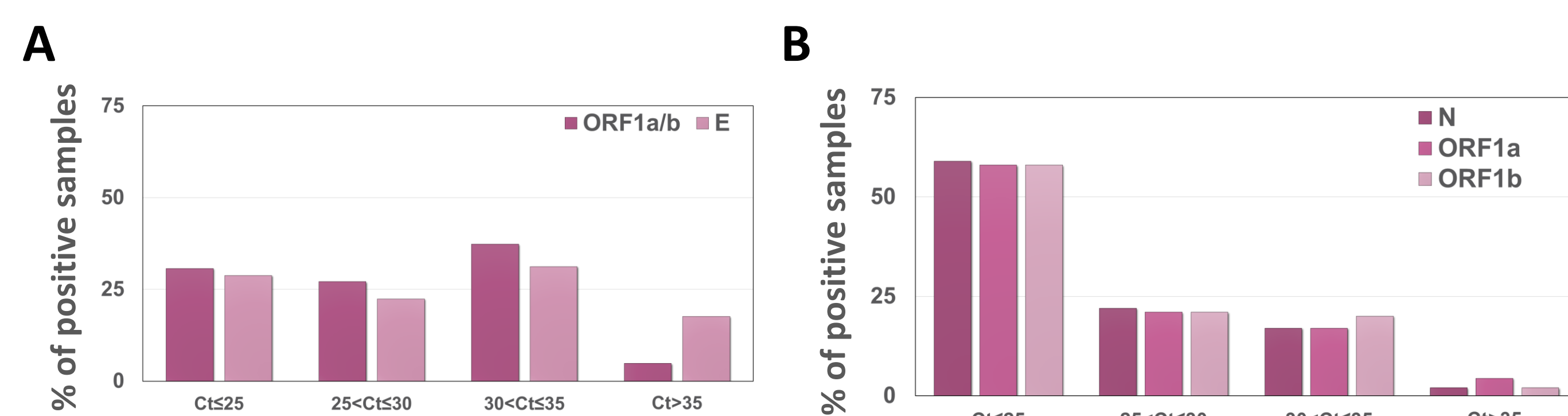


Figure 2. Distribution of Ct values for positive cohort based on A) cobas® SARS-CoV-2 assay and B) TaqPath™ COVID-19 RNase P Combo Kit 2.0\* testing

Of 675 samples, 8 samples were excluded from the analysis due to inconclusive or invalid results. The positive cohort spanned the dynamic range of the assay with a minimum of 20% of samples in each category:  $Ct \leq 25$ ,  $25 < Ct \leq 30$ ,  $30 < Ct \leq 35$  and  $Ct > 35$  (Figure 2). The results showed an excellent concordance between the TaqPath™ COVID-19 RNase P Combo Kit 2.0\* and the cobas® SARS-CoV-2 Assay, with PPA of 97.7% and NPA of 98.2% (Table 1). All 116 VOC samples (Alpha (N=104) and Delta (N=12)) showed a positive result using the TaqPath™ COVID-19 RNase P Combo Kit 2.0 without any inconclusive calls.

Table 1. Concordance between TaqPath™ COVID-19 RNase P Combo Kit 2.0\* and cobas® SARS-CoV-2 Assay

|  |          | cobas® SARS-CoV-2 assay |                  |       |
|--|----------|-------------------------|------------------|-------|
|  |          | Positive                | Negative         | Total |
| TaqPath™ COVID-19 RNase P Combo Kit 2.0* | Positive | 173                     | 9                | 182   |
|  | Negative | 4                       | 481              | 485   |
| Total                                    |          | 177                     | 490              | 667   |
| Positive Percent Agreement (95% CI)      |          | 97.7%                   | (94.3% to 99.4%) |       |
| Negative Percent Agreement (95% CI)      |          | 98.2%                   | (96.5% to 99.2%) |       |

Table 2. Discordant sample resolution using TaqPath™ COVID-19 RT-PCR CE-IVD kit\*

| Sample number | TaqPath™ COVID-19 RNase P Combo Kit 2.0* |       |       |         | Result   | cobas® SARS-CoV-2 Assay | TaqPath™ COVID-19 RT-PCR CE-IVD kit* |
|---------------|--|-------|-------|---------|----------|-------------------------|--------------------------------------|
|               | N  | ORF1a | ORF1b | RNAse P |          |                         |                                      |
| 1             | 31.21                                    | 31.72 | 32.62 | 24.40   | Positive | Negative                | Positive                             |
| 2             | 34.94                                    | 33.74 | 33.84 | 24.81   | Positive | Negative                | Negative                             |
| 3             | 29.34                                    | 29.89 | 30.40 | 23.41   | Positive | Negative                | Negative                             |
| 4             | 31.98                                    | 37.09 | 34.73 | 25.53   | Positive | Negative                | Positive                             |
| 5             | 31.24                                    | 36.41 |       | 20.90   | Positive | Negative                | Positive                             |
| 6             | 28.62                                    | 30.82 | 30.93 | 23.63   | Positive | Negative                | Positive                             |
| 7             | 35.35                                    | 36.65 |       | 22.20   | Positive | Negative                | Negative                             |
| 8             | 27.50                                    | 28.46 | 29.02 | 21.17   | Positive | Negative                | Positive                             |
| 9             | 32.64                                    | 36.17 | 38.22 | 27.97   | Positive | Negative                | Negative                             |
| 10            |  |       |       | 25.48   | Negative | Positive                | Positive                             |
| 11            |  |       |       | 28.71   | Negative | Positive                | Negative                             |
| 12            |  |       |       | 24.49   | Negative | Positive                | Negative                             |
| 13            |  |       |       | 23.73   | Negative | Positive                | Negative                             |

Upon arbitration testing of the 13 discordant samples (Table 2), the sensitivity and specificity of the TaqPath™ COVID-19 RNase P Combo Kit 2.0\* was 99.4% and 99.2% respectively (Table 3).

Table 3. Clinical Sensitivity and Specificity of the TaqPath™ COVID-19 RNase P Combo Kit 2.0\* after discordant sample resolution

|                               |       | TaqPath™ COVID-19 RNase P Combo Kit 2.0 |  |
|-------------------------------|-------|---|--|
| Clinical Sensitivity (95% CI) | 99.4% | (96.9% to 99.9%)                        |  |
| Clinical Specificity (95% CI) | 99.2% | (97.9% to 99.7%)                        |  |

## Conclusion

The highly mutation-resilient TaqPath™ COVID-19 RNase P Combo Kit 2.0\* has excellent concordance with the cobas® SARS-CoV-2 Assay and is a highly accurate method for detection of SARS-CoV-2 including VOC.