

eurogentec.com

Takyon[®] Ultra One-Step **RT-qPCR MM**x For RNA and DNA virus detection



Designed for optimal performance

Further building on its innovative Takyon® line, Eurogentec developed Takyon® Ultra - a unique reverse transcription **qPCR MasterMix** with increased multiplex capabilities and stability. This RT-qPCR kit has been optimized for the sensitive and specific detection of viral RNA and DNA templates such as that from SARS-CoV-2, influenza, Mpox and herpes virus.



RT-qPCR kit Components

- 4X Reaction MasterMix with HotStart Takyon[®] fusion DNA polymerase less sensitive to inhibitors; salts ; dNTPs and stabilizers
- EuroScript II RT PLUS combined with RNAse inhibitor and stabilizers for improved stability and robustness
- ROX™ normalizing dye.

The Advantages of Takyon[®] Ultra

() One-Step RT-qPCR From RNA templates to results in just 1 reaction setup



Multiplex Improved

Co-detect multiple pathogens in a single clinical sample



4X FAST MasterMix

A more concentrated Mix to maximize your template volume, with FAST results delivery



For all Thermocyclers

Individual ROX™ vial included for optimal results on all ROX-based platforms



Low copy detection

Clean negative results and high sensitivity on viral targets.



Automation friendly

24h stability at 20°C for convenient high throughput settings

Proven detection of COVID-19



Highly sensitive viral RNA detection with a broad dynamic range and clean NTCs.

Linear dynamic range in 10-fold dilutions from 500 000 down to 5 copies of SARS-COV-2 RNA Control (TWIST Bioscience #MT007544.1). One-Step Takyon® Ultra Probe 4X MasterMix (yellow) shows a better linear dynamic range than competitor T (dark blue) and better signal than competitor B (light blue) reagents. The amplifications were performed in presence of Yeast T-RNA, under the recommended fast cycling conditions. (NTC: dotted lines)



Takyon® Ultra Performance matches or outperforms most available kits on the market.

Our Optimized Reagents ensure more specific results with clean negative COVID-19 sample results!

Sensitivity and specificity of Sarbeco E gene (SARS-Cov-2 RNA viral target) detection in samples from known COVID-19 positive (patients 1 & 2) or negative (patient 3) patients. The One-Step Takyon® Ultra Probe 4X MasterMix (yellow) shows increased sensitivity compared to Competitor T (dark blue) and Competitor B (light blue) reagents, and increased specificity compared to competitor T, under their recommended fast cycling conditions.



Each component is stable at 20°C for up to 24h providing a convenient solution for automated high-throughput assays set-up.

10-fold linear range from 100 000 down to 10 copies of SARS-COV-2 RNA Control (TWIST Bioscience #MT007544.1) under different storage temperatures for 24h. The One-Step Takyon Ultra Probe 4X MasterMix maintains consistent amplification results after storage for 24h at -20°C (blue, recommended storage conditions), 4°C (grey) or 20°C (yellow) in fast cycling conditions. (NTC: dotted lines)

Takyon® is a registered trademark of Kaneka Eurogentec S.A. ROX™ is a trademark of Applera Corp.

Ordering Information

Cat #	Pack size*	MasterMix Volume
UF-UPRT-C0201	200 – 20 µL RXNs	1 mL
UF-UPRT-C1001	1000 – 20 µL RXNs	5 x 1 mL
UF-UPRT-C1005	5000 – 20 µL RXNs	25 x 1 mL
UF-UPRT-C1010	10,000 – 20 µL RXNs	50 x 1 mL

*Custom batches and packaging available on request

Order on www.eurogentec.com Do you need reagents for commercial use?

Contact us at

custom-qpcr@eurogentec.com