

OPTIMIZED qPCR MASTER MIXES

# Takyon<sup>®</sup> Etheris

Recommended for  
microbial DNA detection

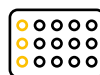
Takyon<sup>®</sup> Etheris is a new generation of qPCR Master Mixes optimized for microbial DNA detection. It combines high sensitivity and robustness with an ultra-purified formulation containing minimal residual host DNA, the main cause of false positives in microbial assays.



**Ultra-low residual  
host DNA**



**Assembled in  
ISO-classified  
cleanrooms**



**Available for SYBR<sup>®</sup>  
& Probe assays**



**Compatible with all  
thermocyclers**

## Optimized for both Probe and SYBR® Green qPCR

Takyon® Etheris is available in Probe and SYBR® Green versions, giving you flexibility for microbial DNA detection workflows.

### Common advantages

- | Compatibility with all qPCR instruments
- | Hot Start Takyon® Fusion DNA polymerase
- | High fluorescence signal

### Additional advantages for probe MMx

- | Enhanced PCR inhibitor tolerance
- | Optimized formulation for high specificity and sensitivity
- | Ideal for multiplexing

## Assembled in cleanroom for trusted qPCR results

Takyon® Etheris is assembled in ISO-classified cleanrooms using ultra-purified reagents with minimal residual host DNA to generate clean No Template Controls (NTC).

Each batch undergoes strict quality control, including background DNA testing, to ensure:

- | Low bacterial nucleic acid content
- | Consistent lot-to-lot performance
- | Reliable results in sensitive qPCR assays



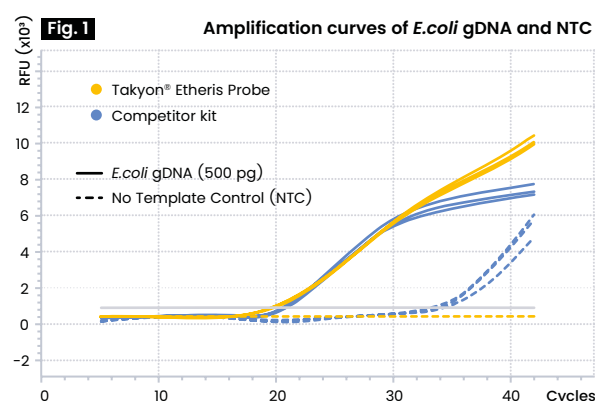


DISCOVER THE  
FULL RESULTS ON  
OUR WEBSITE

## Takyon® Etheris Probe removes background signals

Takyon® Etheris delivers higher sensitivity and stronger amplification signals, with no detectable signal from residual bacterial gDNA. By contrast, the competitor kit produced a Cq signal in the NTC, demonstrating contamination from residual host DNA.

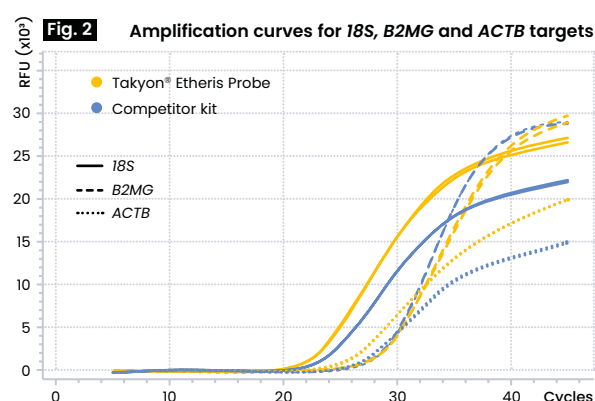
**Fig 1.** Amplification curves with (line) or without (dotted line) *E.coli* gDNA23S (500 pg) in a qPCR reaction using either Takyon® Etheris Probe (orange) or a competitor kit (blue). Each assay was performed in technical triplicates. All reactions were run according to the respective manufacturer's protocols.



## Outperforms specialized competitor's reagents

Compared to the competitor kit Takyon® Etheris demonstrates improved sensitivity and stronger fluorescence signal intensity across all targets, proving its value in multiplex conditions.

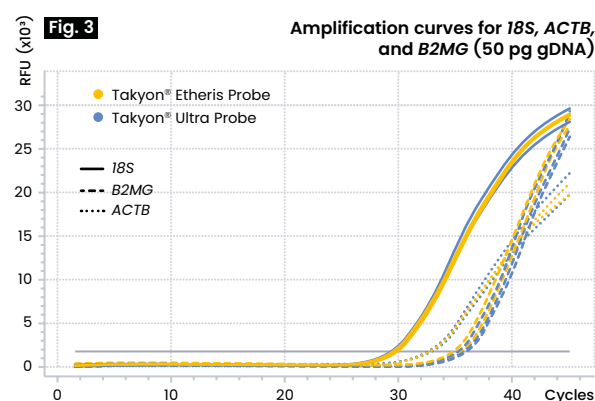
**Fig 2.** Amplification curves obtained with Takyon® Etheris Probe qPCR Master Mix (orange) and a competitor kit (blue) in a triplex assay targeting *18S*, *B2MG*, and *ACTB* genes. Each reaction was performed in triplicate using 5 ng of DNA template according to the respective manufacturer's protocols.



## Keeps the performance of the Takyon® range

Takyon® Etheris and Takyon® Ultra exhibit similar sensitivity and amplification performance across all three tested targets at 50 pg of human gDNA (*18S*, *ACTB* and *B2MG*). Mean Cq values are closely aligned between the two formulations, confirming their equivalent efficiency in detecting low-copy templates.

**Fig 3.** Amplification curves obtained with Takyon® Etheris Probe qPCR Master Mix (orange) and Takyon® Ultra Probe qPCR Master Mix (blue) for amplification of three human targets: *18S*, *ACTB*, and *B2MG*. Each reaction was performed in triplicate using 50 pg of human genomic DNA as template.



## For any assay aimed at detecting microbial DNA



### Environmental microbiology

Reliable results even in complex soil or plant samples.



### Water quality monitoring

Detect microbes in drinking or wastewater with low background.



### Food safety & quality

Ensure pathogen-free products and validated hygiene control.



### Clinical & diagnostic microbiology

Trust your molecular tests with clear, contamination-free amplification.



### Microbiome analysis

Profile microbial diversity with confidence and precision.



### DNA barcoding

Track and identify species across your production chain.

## What will you find in your kits?

	Takyon® Etheris for SYBR®-based assays	Takyon® Etheris for Probe-based assays
<b>2X Master Mix</b> with Hot Start Takyon® fusion DNA Polymerase	Blue Master Mix	Clear Master Mix
<b>ROX Normalizing dye</b> provided separately making the kit compatible with all thermocyclers	1 tube of 0.5 mL	1 tube of 0.5 mL
<b>Available Size</b> number of 20 µL reactions	150 RXNs – 750 RXNs 5 packs of 750 RXNs 10 packs of 750 RXNs	150 RXNs – 750 RXNs 5 packs of 750 RXNs 10 packs of 750 RXNs



SCAN TO ORDER  
A FREE SAMPLE



SCAN TO ORDER  
A FREE SAMPLE

Do you need larger reagent volumes  
for routine or commercial use?

Scan this code and contact us at  
[custom-qpcr@eurogentec.com](mailto:custom-qpcr@eurogentec.com)

