

Takyon® Etheris

Probe 2X qPCR Master Mix^[1]

UF-EPMT-C0101

[1.5 mL]
[150 x 20 µL RXNs]

UF-EPMT-C0701

[7.5 mL]
[750 x 20 µL RXNs]

UF-EPMT-C0705

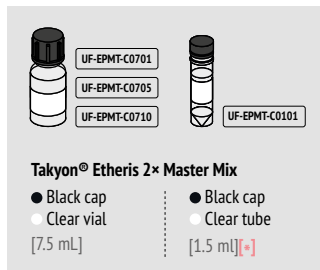
[5 packs of 7.5 mL]

UF-EPMT-C0710

[10 packs of 7.5 mL]

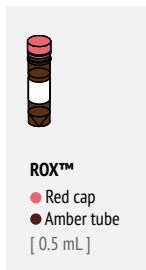
Kit content (for 750 (150^[*]) - 20 µL reactions)

INCLUDED



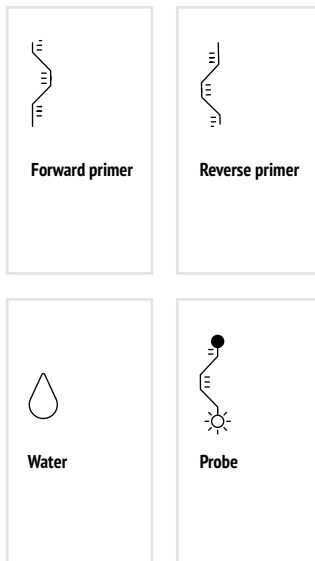
- Takyon® HotStart DNA polymerase
- MgCl₂ (5.5 mM final concentration)
- dNTPs
- Stabilizers

OPTIONAL



For signal
normalization

NOT INCLUDED



Storage

Takyon® kit must be stored in a constant temperature freezer in the dark.

SHORT TERM STORAGE

⌚ **6** Month stability

🌡️ **4°C**

LONG TERM STORAGE

⌚ **24** Month stability

🌡️ **-15°C**
-25°C

Optimization tips

Refer to the primer and probe design guidelines, custom assay design recommendations, primer titration matrix, and primer-probe ratio matrix for best results.

For new users of this kit, conducting a primer titration matrix may be required to ensure optimal performance.



**TECHNICAL
INFORMATION**

[1] Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

⚠ ROX™ Adjustment before first use

IMPORTANT

Before first use, convert your Master Mix to the correct ROX™ level.

Table 1 shows how much ROX™ passive reference to add to Takyon® Etheris qPCR Master Mix, according to the thermocycler used.

No ROX™ is needed for most thermocyclers not listed.

| qPCR PLATFORM | 1.5 ML MMX [ROX™ TO ADD] | 7.5 ML MMX [ROX™ TO ADD] |
|---|-----------------------------|-----------------------------|
| Mx3000P®, 3005P, 4000 | 12 µL | 60 µL |
| ABI Prism® 7500, FAST 7500, ViiA7™, QuantStudio™ | 4.2 µL | 21 µL |
| ABI Prism® 5700, 7000, 7300, 7700, 7900, FAST 7900, ABI Step One, One Plus MasterCycler® ep realplex I & II (rev. =< 2.1) | 60 µL | 300 µL |

Table 1 : ROX™ addition according to qPCR Platform.

Recommended protocol

1 Thaw all required reagents completely and put them **on ice**.



2 Prepare the REACTION MIX in excess to correct for dispensing losses (e.g. a 100-reactions mix for 96-reactions). **Add all components together**, except for the template. Mix thoroughly by pipetting or inversion. **Spin down**.



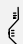


| REACTION MIX COMPONENTS | VOLUME / REACTION | FINAL CONCENTRATION |
|---|-------------------|---|
|  Takyon® MMX | 5 µL | 1× |
|  Forward primer | 2 µL | 50-900 nM[3] |
|  Reverse primer | 2 µL | 50-900 nM[3] |
|  Probe | 2 µL | 100-250 nM[3] |
|  RNase-free Water | x µL[2] | (volume is 20 µL minus all other components)[4] |
| Total Mix/Reaction | 20 µL | |

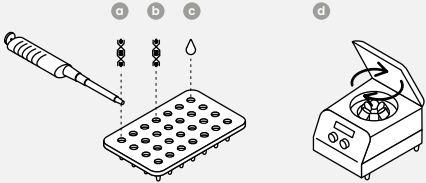
Table 2 : Reaction mix preparation.

3 Add the REACTION MIX to individual reaction vials.



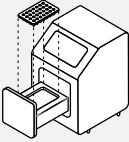
4 Pipette into your qPCR plate/tubes either[4]:

- a 2.5 µL of template cDNA / DNA
- b 2.5 µL of positive control
- c 2.5 µL of water / buffer for negative control



d Gently mix on a magnetic stirrer and centrifuge to avoid bubbles.

5 Program the Real-Time thermocycler using the following recommended parameters:



Takyon® Etheris produces consistent and sensitive results under FAST[5] cycling conditions.

| | °C | TIME |
|---------------------------------|------|-------------|
| Takyon® activation | 95°C | 3 min |
| 40 CYCLES[5] | | |
| Denaturation[6-7] | 95°C | 15 sec |
| Annealing / extension[7] | 60°C | 20 - 30 sec |

Table 3 : Cycling conditions.

Trademarks : ABI Prism® 5700, 7000, 7500, 7700 and 7900 are registered trademarks of Applied Biosystems, Inc. | MasterCycler® is a registered trademark of Eppendorf, Inc. | Mx3000P®, Mx3005P® and Mx4000® are registered trademark of Stratagene | ROX™ is a trademark of Applied Biosystems. | Takyon® is a registered trademark of Kaneka Eurogentec S.A.

[2] Adjust water so total volume is 20 µL after adding all components. [3] Start with primers at 300 nM and probe at 250 nM; optimize if needed. [4] Mix 17.5 µL reaction mix with 2.5 µL DNA for a 20 µL final volume; adjust if required. [5] Use FAST cycling only on FAST-block cyclers; prefer amplicons <120 bp. [6] For complex templates (e.g., plant/genomic DNA), increase denaturation time. [7] Set annealing temperature according to primer Tm. For longer/difficult templates, extend annealing/extension to 40 s. Some FAST cyclers allow shorter denaturation/annealing for quicker qPCR; if amplification is poor, lengthen extension in 5 s steps.