

Takyon® Etheris SYBR® 2X qPCR Master Mix Blue^[1]

UF-ESMT-B0101

[1.5 mL]
[150 x 20 µL RXNs]

UF-ESMT-B0701

[7.5 mL]
[750 x 20 µL RXNs]

UF-ESMT-B0705

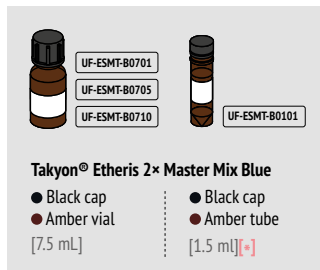
[5 packs of 7.5 mL]

UF-ESMT-B0710

[10 packs of 7.5 mL]

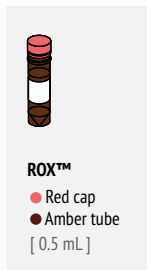
Kit content (for 750 (150^[*]) - 20 µL reactions)

INCLUDED

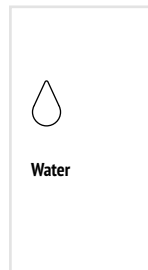
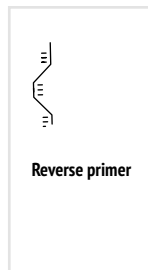
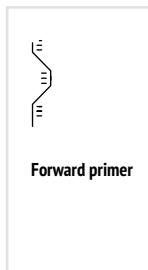


- Takyon® HotStart DNA polymerase
- SYBR® Green
- Inert Blue Dye
- MgCl₂ (2.5 mM final concentration)
- dNTPs
- Stabilizers

OPTIONAL

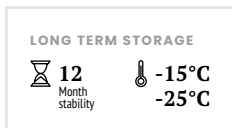
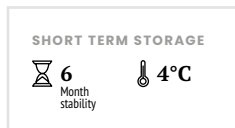

For signal
normalization

NOT INCLUDED



Storage

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



Optimization tips

Refer to the primer design guidelines, custom assay design recommendations and primer titration 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



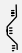

Table 1 : ROX™ addition according to qPCR Platform.

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	50-300 nM[3]
 Reverse primer	2 µL	50-300 nM[3]
 RNase-free Water	x µL[2]	(volume is 20 µL minus all other components)[4]
Total Mix/Reaction	20 µL	


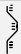


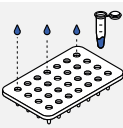
REACTION MIX COMPONENTS	VOLUME / REACTION	FINAL CONCENTRATION
 Takyon® MMx	5 µL	1×
 Forward primer	2 µL	50-300 nM[3]
 Reverse primer	2 µL	50-300 nM[3]
 RNase-free Water	x µL[2]	(volume is 20 µL minus all other components)[4]
Total Mix/Reaction	20 µL	

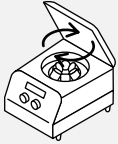
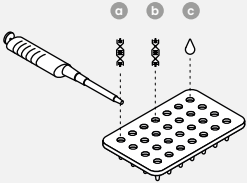
Table 2 : Reaction mix preparation.

3 Add the REACTION MIX to individual reaction vials.



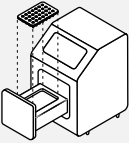
4 Pipette into your qPCR plate/tubes either[4]:

- 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.

	T°C	TIME
Takyon® activation	95°C	3 min
40 CYCLES[5]		
Denaturation[6]	95°C	3 - 10 sec
Annealing / extension[7]	60°C	20 - 60 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 Applera Corp. | SYBR® Green is a registered trademark of Life Technologies Corp. | 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 100 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.