

---

 Technical **Data Sheet**


---

## TAKYON® One-Step Kit Converter (Euroscript II + Additive) UF-RTAD-D0701 kit • UF-RTAD-D0101 sample

Eurogentec products are sold for research or laboratory use only and are not to be administered to humans or used for medical diagnostics.

### Kit contents

The standard kit contains enough reagents to convert 7.5 mL of 2X Takyon® qPCR dTTP MasterMix or 3.75 mL of 4X Takyon qPCR MasterMix into a One-Step qPCR kit. It contains:

Description	Volume	Details
	UF-RTAD-D0701 (UF-RTAD-D0101)	
Euroscript II RT / RNase inhibitor (white cap)	150 µL (30µL)	50 u/µL RT 20 u/µL RNase inh.
Additive (blue cap)	150 µL (30µL)	Improve results on some viral and FFPE samples

### Storage conditions

For long term storage the Euroscript II RT should be stored at -65 °C to -75 °C in a constant temperature freezer. When stored under these conditions the RT is stable for 2 years. The Euroscript II RT and additive can be stored between -15°C and -25°C for up to 12 months.

### Intended use

The reagents have been optimized for use in One-Step assays with Takyon® qPCR dTTP MasterMixes.

### Procedure

- 1 Thaw all required reagents, incl. Takyon® qPCR MasterMix, completely and put them on ice, except for the EuroScript II RT and RNase Inhibitor, which should be kept in the freezer until required for use. Mix all reagents well by inversion and spin them down prior to pipetting.
- 2 Prepare the reaction mix. To correct for dispensing losses, prepare an excess of the mix (e.g. a 100-reaction mix for 96 reactions). Reaction set up should be done on ice.

– Add all components together, except for the template, as per Table 2 in the Takyon® MasterMix Technical Data sheet.

- Add Takyon® MasterMix (2X or 4X)
- Add primers & probe(s) at optimized concentration
- Add 0.20µl of Euroscript II RT per 20µl reaction (final volume)
- *Optional: add 0.20µl of additive per 20µl reaction (final volume). Additive improves results on some viral templates and on FFPE samples.*
- Add RNase-free water to reach final volume

– Mix thoroughly by inversion. Spin down.

3 Add the reaction mix to individual reaction vials.

4 Add the template to individual reaction vials, gently mix on a magnetic stirrer and centrifuge to avoid bubbles. Negative control containing no RNA template should always be included. Optionally, a no RT-control should be set up in tubes / wells, which does not contain the EuroScript II RT/RNase Inhibitor.

5 The Takyon® One-Step Converter reagents will produce consistent and sensitive results under FAST and REGULAR cycling conditions. Program the Real-Time thermocycler using the following recommended cycling parameters:

#### **A. Reverse transcription step: 10 min. at 48 °C**

For difficult templates, increase RT step by increment of 10', up to a total of 30', to improve reaction yield.

#### **B. c-DNA amplification step:** please follow the recommended Takyon MasterMix cycling parameters (see table 3 of the Takyon Technical Data Sheet).

**For further information please contact our Customer Help Desk:**

#### For Europe:

E-mail: info@eurogentec.com

Tel: +32 4 372 76 65 • Toll-free: + 800 666 00 123