



Takyon™ dry

Dry MasterMixes for qPCR and RT-qPCR

NEW
product

For Life Science Research

⇒ LYOPHILISED FORMAT

Takyon™ Dry is a new lyophilised ready-to-use MasterMix format for real-time qPCR and RT-qPCR Probe assays.

The dried reagent contains the new robust **Takyon™ fusion enzyme**. Takyon™ Dry produces consistent and sensitive results under regular and fast cycling protocols.

Takyon™ Dry reagents are compatible with all qPCR platforms in a ready-to-use format (ROX, low ROX and no ROX). ■

*Ship it, store it,
use it anywhere,
anytime*



BENEFITS



Eco-friendly & Economical

Takyon™ Dry reagents are environmental friendly. Ambient shipping allows for lightweight, dry ice- and polystyrene foam-free parcels while room temperature storage circumvents the need for a refrigerator.



Stable at room temperature

The lyophilisation process makes the reagents stable for 18 months at ambient temperature, greatly simplifying storage and handling of the kits.



Efficient

Make no compromise on performance! Takyon™ Dry reagents demonstrate equivalent performance to that of the efficient and sensitive in-solution DNA and RNA Takyon™ MasterMixes.



Flexible

Takyon™ Dry mix can be easily and instantly reconstituted in any volume up to a 5X concentrated MasterMix.

Takyon™ dry

Takyon™ qPCR MasterMixes for Probe Assays

Takyon™ RT-qPCR MasterMixes for One-Step Probe Assays



ORDERING INFORMATION

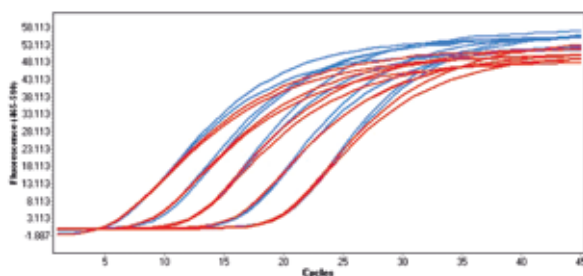
Product	Quantity	Cat.#
Takyon™ Dry Rox Probe MasterMix dTTP	3x50 rxn	UFD-RPMT-C0101
Takyon™ Dry Low Rox Probe MasterMix dTTP	3x50 rxn	UFD-LPMT-C0101
Takyon™ Dry No Rox Probe MasterMix dTTP	3x50 rxn	UFD-NPMT-C0101



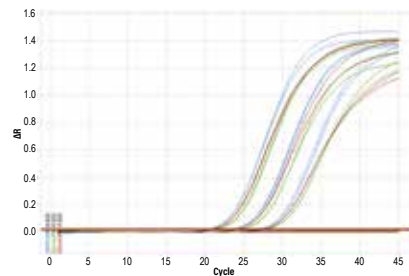
ORDERING INFORMATION

Product	Quantity	Cat.#
Takyon™ Dry Rox One-Step RT Probe MasterMix	3x50 rxn	UFD-RPRT-C0101
Takyon™ Dry Low Rox One-Step RT Probe MasterMix	3x50 rxn	UFD-LPRT-C0101
Takyon™ Dry No Rox One-Step RT Probe MasterMix	3x50 rxn	UFD-NPRT-C0101

Wet and Dry qPCR/ RT-qPCR Kit performance are equivalent
Other formulations are available on request.



Detection of dilution series of 28S rRNA on a LC480 using Takyon™ Dry (red) or catalogue UFN5X-RT0501 Takyon™ 5X One-Step MMx (blue) under identical cycling conditions.



Detection of dilution series of MS2 RNA assay on a ECO cycler using Takyon™ Dry resuspended in buffer S (red), in water (green) or catalogue UFN5X-RT0501 Takyon™ 5X One-Step MMx (blue) under identical cycling conditions.

HOW TO ORDER

WWW.EUROGENTEC.COM



ONLINE

Every Takyon™ MasterMix can be ordered online. Log in to your Eurogentec account and select the products of your choice.

EMAIL ORDERING

Please send your request to order@eurogentec.com

QUOTATION

For more information or to get a quotation please contact your local sales representative
<http://www.eurogentec.com/sales-contact.html>
or isales@eurogentec.com

www.eurogentec.com/qpcr-takyon-dry.html

CUSTOM FORMULATION

Custom specification and custom kit lyophilisation can be discussed, please contact us at:
custom-pcr@eurogentec.com