

EUROPE

KANEKA EUROGENTEC SA - LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00 • Fax: +32 4 372 75 00 Toll-free: + 800 666 00 123 • info@eurogentec.com • www.eurogentec.com

NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560 Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • service@anaspec.com • www.anaspec.com



Experience true partnership

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

Technical Data Sheet

Takyon™ BLUE Additive

UF-BLUE-ADD

[3750 (20 µL) RXN]

Description

The Takyon™ BLUE Additive is a non-fluorescent inert dye which considerably enhances the contrast between reagent and plasticware, making verification of MasterMix dispensing quick, easy and foolproof.

Storage condition

For long term storage the Takyon™ BLUE Additive should be stored at -20 °C in a constant temperature freezer. When stored under these conditions the reagent is stable for 2 years upon batch release.

The Takyon™ BLUE Additive can be stored at 4°C for up to 3 months.

Avoid repetitive freeze-thaw cycles.

Product content

The Takyon™ BLUE Additive contains enough reagents to convert up to 3750 - 20 µL Takyon™ qPCR reactions.

Reagent	Volume
Takyon™ BLUE Additive (Blue cap)	1 x 375 µL

Recommendations of use

Convert Takyon™ Mastermixes upon first use as per follow:

1. Thaw the Takyon™ MasterMix and the Takyon™ BLUE additive
2. Add the Takyon™ BLUE additive to the Takyon™ MasterMix vial as per follow:
 - 15 µL into 1.5 mL of 2X MasterMix
 - 15 µL into 0.6 mL of 5X MasterMix
 - 25 µL per 1 mL tube of 5X MasterMix
 - 75 µL per 7.5 mL bottle of 2X MasterMix
3. Gently homogenise the Takyon™ MasterMix by swirling the vial
4. Follow MasterMix storage recommendation for the converted kit

Takyon™ is a trademark of Kaneka Eurogentec

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

For USA:

E-mail: service@anaspec.com

Tel.: +1-510-791-9560 • Toll-free: +1-800-452-5530