

EUROPE

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

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

Uracil-N-Glycosylase (UNG)/dUTP Mix

RT-UTPUNG-005 • RT-UTPUNG-020 • RT-UTPUNG-033

Storage conditions

For long term storage the dUTP/UNG blend should be stored at -20°C in a constant temperature freezer. When stored under these conditions the reagent is stable for 2 years. Avoid repetitive freeze-thaw cycles.

Source

Recombinant *E.coli* UNG purified from an *E.coli* expressing clone.

Description

E.coli Uracil-N glycosylase (UNG) hydrolyses the release of free uracil from uracil-containing DNA. UNG efficiently hydrolyses single or double-stranded DNA, but not from oligomers (6 or less bases).

Table 1

| Reagent | Reference | Volume |
|----------------|---------------|--------|
| dUTP/UNG Blend | RT-UTPUNG-005 | 40 µL |
| | RT-UTPUNG-020 | 200 µL |
| | RT-UTPUNG-033 | 335 µL |

Procedure

Convert a dTTP MasterMix kit before first use:

- Thaw the MasterMix and the vial containing dUTP/UNG at RT
- Add the dUTP/UNG blend to the MasterMix vial as follow:
- 40 µL into 1.5 mL of 2X MasterMix
- 40 µL into 0.6 mL of 5X MasterMix
- 67 µL per 1 mL tube of 5X MasterMix
- 200 µL per 7.5 mL bottle of 2X MasterMix
- Gently homogenize the MasterMix by swirling the vial
- Follow MasterMix storage recommendation for the converted kit

Convert a qPCR reaction into a carryover ready assay (for Core kit):

- Thaw all required reagents
- Add 0.25 µL of dUTP/UNG per 20 µL reaction (final volume - refer to the TDS of the corresponding Core kit)
- Gently homogenize the Mix by pipetting or inverting the tube a few times. Spin down.
- Start the reaction

For further information please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com

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

For USA:

E-mail: service@anaspec.com

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