

Uracil-N-Glycosylase (UNG)/dUTP Mix

• RT-UTPUNG-020 •

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 (µL)
dUTP/UNG Blend	RT-UTPUNG-020	200

Procedure

It is essential to avoid using UNG in PCR cycles where the annealing/extension temperatures are below 55°C to ensure optimal and consistent results. Temperatures of at least 55°C should be used throughout the cycling protocol to avoid degrading the PCR products.

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
 - 56 µL per 1 mL tube of 4X 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

Further information available through Eurogentec web site : www.eurogentec.com.

For further information please contact our Customer Help Desk:

E-mail: info@eurogentec.com
Tel: +32 4 372 76 65

FOR RESEARCH USE ONLY NOTICE TO PURCHASER: LIMITED LICENSE
The sale of this product is conditioned on the buyer using the product only in research conducted by the buyer, and not to sell or otherwise commercialize this product or its components (1) in manufacturing; (2) to provide a service, information, or data in exchange for payment; (3) for therapeutic, diagnostic or prophylactic purposes; or (4) for resale, not even if resold for use in research. For information on purchasing a license to this product for purposes other than research, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, CA 92008 USA or outlicensing@lifetech.com.