

Uracil-DNA Glycosylase Technical Data Sheet

Reference: RT-0610-03 / RT-0610-15

Products and procedures describe in this protocol are intended for research purposes only. Not for use in diagnostic procedures for clinical purposes. **FOR IN VITRO USE ONLY**

Storage conditions

For long term storage the Passive reference should be stored at -15 °C to -25 °C in a constant temperature freezer. When stored under these conditions the reagents are stable for 2 years. Avoid repetitive freeze-thaw cycles.

Source

Purified from *E.coli* expressing the *ung* gene of *E.coli* on a plasmid.

Description

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

Concentration

RT-0610-03: 1,000 units per ml
RT-0610-15: 5,000 units per ml

Supplied with a vial of 10x reaction buffer (200mM Tris-HCl, (pH8.0), 10mM EDTA, 10mM DTT)

Contents

RT-0610-03: 300 U @ 1u/μl
RT-0610-15: 1,500 U @ 1u/μl

Unit definition

One unit is defined as the amount of enzyme that catalyses the release of 60pmol of uracil per minute from double-stranded, uracil containing DNA.

Procedure when using any PCR or qPCR kits containing dUTP

Uracil-N-Glycosylase can be added to a final concentration of 0.01 u/μl:
0.5 μl of 1 u/μl UNG per 50 μl reaction

Protocol

If UNG is required, please add an initial step before the usual cycling protocol: 2 min at 50 °C.

(Hold at 50 °C when the cycling protocol is finished)