

Eurogentec Headquarters

LIEGE SCIENCE PARK • 4102 SERAING • Belgium • Tel: +32 4 372 76 65
Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com

Eurogentec North America

3347 Industrial Ct • Suite A • San Diego, CA 92121 • USA • Tel: +1 858 793 26 61 Fax: +1 858 793 26 66 • info.usa@eurogentec.com • www.eurogentec.com

Fluorescein Additive Technical Data Sheet

Reference: RT-FLUO-ADD

Products and procedures describe in this protocol are intended for research purposes only.

Storage conditions

For long term storage the Fluorescein Additive should be stored at -15 °C to -25 °C in a constant temperature freezer. When stored under these conditions the reagents are stable for a minimum of twelve months. Avoid repetitive freeze-Thaw cycles.

Contents

The Fluorescein Additive contains 1 ml of 1mM fluorescein in DMSO.

Procedure

To enable the use of the qPCR Core kit for SYBR® green I No ROX/RT-SN10-05NR or the qPCR MasterMix Plus for SYBR® green I No ROX / RT-SN2X-03+NR on an iCycler iQ® or My iQ, the addition of fluorescein is required.

- Prepare a working solution by adding 1 μ I of fluorescein additive to 999 μ I PCR grade water (1:1000 dilution). This working solution can be aliquoted and stored at -20 °C. Avoid Freeze-thaw cycles.
- Dilute this working solution 1:100 in the final reaction mix, so for example add 0.5 µl of the working solution per 50 µl reaction.