

PRODUCT INFORMATION

ROX Solution, 50 μ M

#R1371 1 x 1 mL

Lot _ Expiry Date _

Store at -20°C

www.thermoscientific.com/onebio

Description

ROX Solution contains ROX dye as a 50 μ M aqueous solution at pH 7.6.

The excitation maximum and emission maximum of ROX are at 580 nm and 621 nm, respectively.

Applications:

For use in qPCR as an internal reference when normalizing the fluorescent signal with instruments which can detect ROX. For detailed recommendations refer to the manual of qPCR.

CERTIFICATE OF ANALYSIS

Functionally tested in real-time PCR in parallel 20 μ L reactions containing 10-fold dilutions of human genomic DNA to demonstrate linear resolution over five orders of dynamic range.

Quality authorized by:  Jurgita Zilinskiene

NOTICE TO PURCHASER: DISCLAIMER OF LICENSE

No license is conveyed with the purchase of this product under US Patent No. US 5,928,907 and corresponding claims in patents outside the United States relating to the use of a fluorescent indicator as a passive reference in real time PCR. For further information contact the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals.

Please refer to www.thermoscientific.com/onebio for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Rev.1