

PRODUCT INFORMATION

40X Yellow Sample Buffer

#R1381 4 × 1.25 mL

Lot _

Store at -20°C

Description

40X Yellow Sample Buffer is an aqueous solution of the inert yellow dye at pH 8.0.

The absorption maximum of the yellow dye is at 413 nm.

Applications

For use with Luminaris Color qPCR Master Mixes to color and track the samples. For detailed recommendations refer to the manual of qPCR.

CERTIFICATE OF ANALYSIS

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

Quality authorized by:



Jurgita Zilinskiene

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