



ROX Reference Dye

Reference dye for fluorescence signal normalization

Cat. No.	Amount
PCR-351	1 ml (25 µM)

For *in vitro* use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: store dark

Shelf Life: 12 months

Form: liquid, supplied in 20 mM Tris-HCl pH 8.5, 0.1 mM EDTA and 0.01 % Tween-20.

Color: pink

Concentration: 25 µM (50x conc.)

Spectroscopic Properties: λ_{exc} 576 nm, λ_{em} 601 nm

Description:

ROX reference dye is optimized for block-based real-time PCR instruments that are compatible with the evaluation of the ROX reference signal. The dye does not take part in the PCR reaction but allows to normalize for non-PCR related signal variation and provides a baseline in multiplex reactions. The dye is supplied as 50x concentration.

Preparation of the qPCR master mix:

Preparation of a master mix is crucial in quantitative PCR reactions to reduce pipetting errors. Vortex the ROX reference dye thoroughly before use. Add ROX reference dye to your PCR assay as indicated below.

final ROX concentration	20 µl PCR assay	50 µl PCR assay
low ROX assay 50 nM	0.04 µl	0.1 µl
high ROX assay 500 nM	0.4 µl	1.0 µl

Related Products:

qPCR Core Kits
Dual labeled fluorescent probes
Custom primers