



## ROX Reference Dye - 100 $\mu$ M

Reference dye for fluorescence signal normalization

Cat. No.	Amount
PCR-356-1ML	1 ml (100 $\mu$ M)
PCR-356-10ML	10 ml (100 $\mu$ M)
PCR-356-100ML	100 ml (100 $\mu$ M)

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** store dark

**Shelf Life:** 12 months

**Form:** liquid

**Color:** pink

**Concentration:** 100  $\mu$ M

**Spectroscopic Properties:**  $\lambda_{exc}$  576 nm,  $\lambda_{em}$  601 nm

### Description:

ROX reference dye is recommended for block-based real-time PCR instruments that are compatible with the evaluation of a 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 100  $\mu$ M concentration.

### Content:

100  $\mu$ M ROX dye in 20 mM Tris-HCl pH 8.5, 0.1 mM EDTA and 0.01 % Tween-20

### Preparation of the qPCR master mix with ROX

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

assay set-up	final ROX concentration
low ROX assay	50 nM
high ROX assay	500 nM

### Related Products:

qPCR Core Kits  
Dual labeled fluorescent probes  
Custom primers