Glowing products for science

ROX Reference Dye, 25 uM in TE Buffer

A dye used to normalize well-to-well variations that may occur during quantitative, real-time PCR reactions.



Product attributes

 Storage Conditions
 Store at -10 to -35 °C

 Excitation/Emission
 575/600 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

ROX reference dye is used to normalize well-to-well variations that may occur during quantitative, real-time PCR reactions.

- Red dye used to normalize qPCR reactions
- Won't interfere with PCR reactions
- Supplied as a 25 uM solution

References

1. J Feline Med Surg (2019) Feb;21(2):124-132 doi: 10.1177/1098612X18762017

This datasheet was generated on November 4, 2025 at 07:51:56 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rox-reference-dye-25-um-in-te-buffer/