



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 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

## 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](https://doi.org/10.1177/1098612X18762017)

This datasheet was generated on January 6, 2026 at 09:09:26 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/rox-reference-dye-25-um-in-te-buffer/>