

AMPIGENE® qPCR

Probe Mix with Separate ROX

Optimized probe-based qPCR

AMPIGENE® qPCR Probe Mix with Separate ROX uses the latest developments in polymerase technology and buffer chemistry to enhance qPCR speed, yield, and specificity. AMPIGENE® qPCR Probe Mix with Separate ROX uses advanced hot-start technology for superior sensitivity. The enzyme and buffer system allow for superior qPCR performance on a broad range of templates including complex templates such as mammalian genomic DNA.

AMPIGENE® qPCR Probe Mix with Separate ROX contains AMPIGENE® Hot Start *Taq* DNA Polymerase, a robust enzyme for all your everyday PCR applications, dNTPs, MgCl, DMSO, and propriety enhancers, enabling high sensitivity and high-fidelity qPCR of a wide range of targets and fragment sizes. The buffer system allows the efficient amplification of GC-rich and AT-rich templates under standard cycling conditions. AMPIGENE® qPCR Probe Mix with Separate ROX is engineered for use with a wide range of probe technologies such as TaqMan® and molecular beacon probes.

The AMPIGENE® ROX passive reference dye is supplied in a separate vial at 50X concentration.

Ordering Information

[Order Online »](#)

ENZ-NUC127-1000	1000Reactions
-----------------	---------------

Manuals, SDS & CofA

[View Online »](#)

- Hot start technology prevents primer-dimers formation, improving sensitivity and specificity
- Better performance for amplification of GC-rich targets compared to competitors' product
- Separate vial of ROX allows for flexibility of usage. The qPCR mix can be used without the reference dye ROX, with low concentration of ROX, and for high ROX amplification
- Consistent and reliable lot-to-lot performance, delivering reproduce results

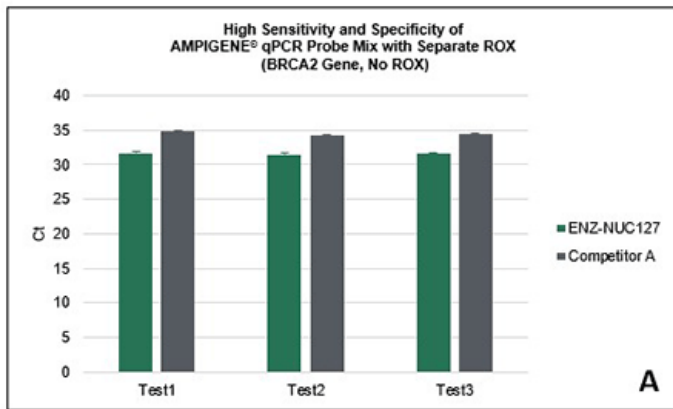


Figure 1A. Amplification results for the breast cancer gene 2 (BRCA2) were obtained from human genomic DNA samples in absence of ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Probe Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

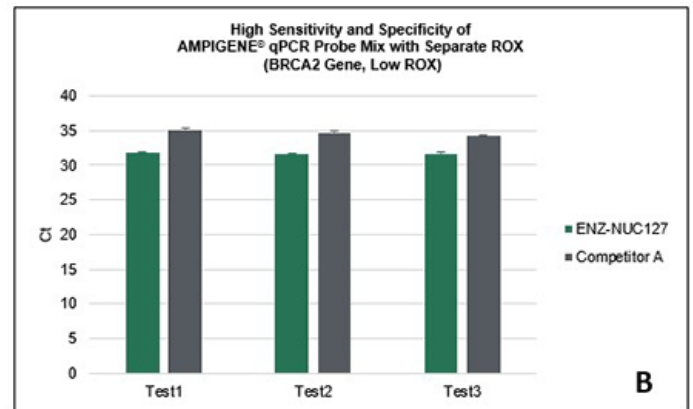


Figure 1B. Amplification results for the breast cancer gene 2 (BRCA2) were obtained from human genomic DNA samples with low ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Probe Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

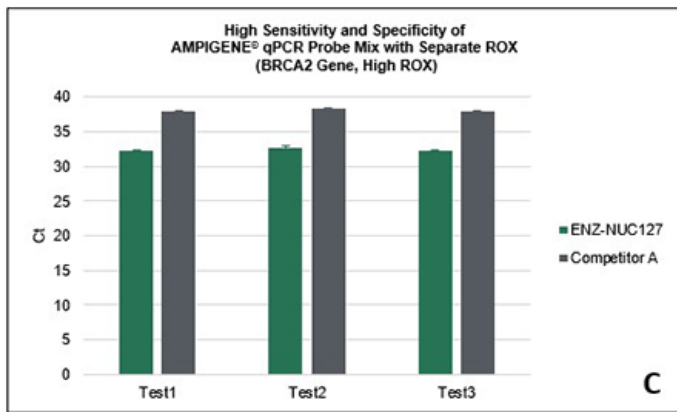


Figure 1C. Amplification results for the breast cancer gene 2 (BRCA2) were obtained from human genomic DNA samples with high ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Probe Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

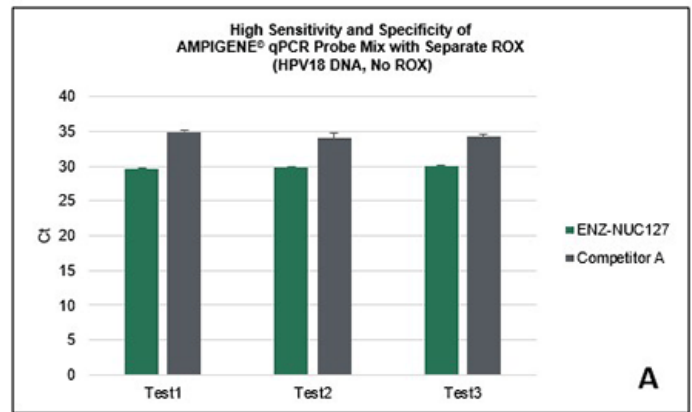


Figure 2A. Amplification results for the human papillomavirus type 18 (HPV18) DNA sample were obtained from human genomic DNA samples in absence of ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Probe Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

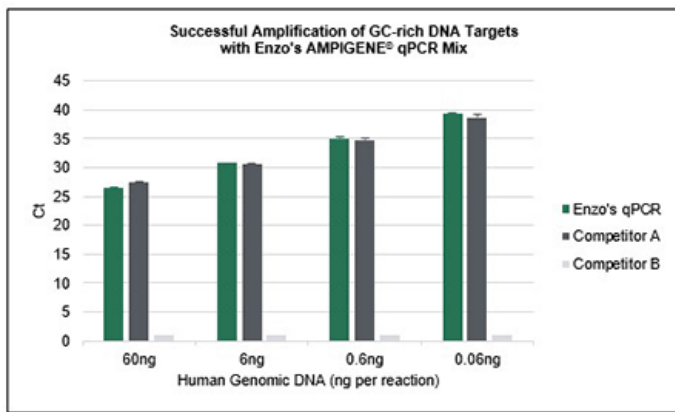


Figure 3. Amplification results for the GC-rich DNA target were obtained by targeting a human genomic region that contains 72.9% GC content. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system. The reactions with competitors' reagents were included and the performances were assessed. Product from competitor B failed the amplification.

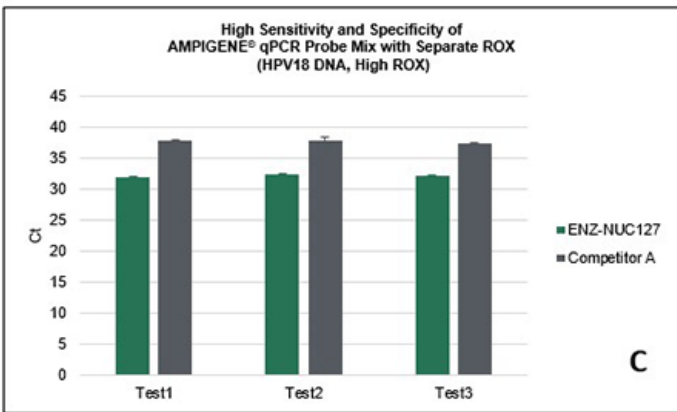


Figure 2C. Amplification results for the human papillomavirus type 18 (HPV18) DNA sample were obtained from human genomic DNA samples with high ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Probe Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

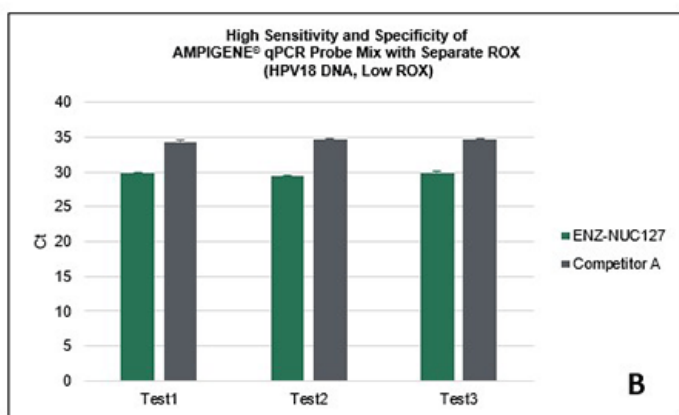


Figure 2B. Amplification results for the human papillomavirus type 18 (HPV18) DNA sample were obtained from human genomic DNA samples with low ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Probe Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application qPCR

Application Notes For probe-based qPCR

Contents 2X AMPIGENE® qPCR Probe Mix with Separate ROX
(AMPIGENE® Hot Start *Taq* DNA Polymerase, MgCl₂,
DMSO, dNTPs, and enhancers)
50X AMPIGENE® ROX

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)