

AMPIGENE® qPCR

Green Mix with Separate ROX

Optimized dye-based qPCR

AMPIGENE® qPCR Green Mix with Separate ROX uses the latest developments in polymerase technology and buffer chemistry to enhance qPCR speed, yield, and specificity. AMPIGENE® qPCR Green Mix with Separate ROX uses advanced hot-start technology for superior sensitivity. AMPIGENE® qPCR Green Mix with Separate ROX contains AMPIGENE® Hot Start *Taq* DNA Polymerase enzyme, PCR green dye, 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. The AMPIGENE® ROX passive reference dye is supplied in a separate vial at 50X concentration.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-NUC126-1000	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

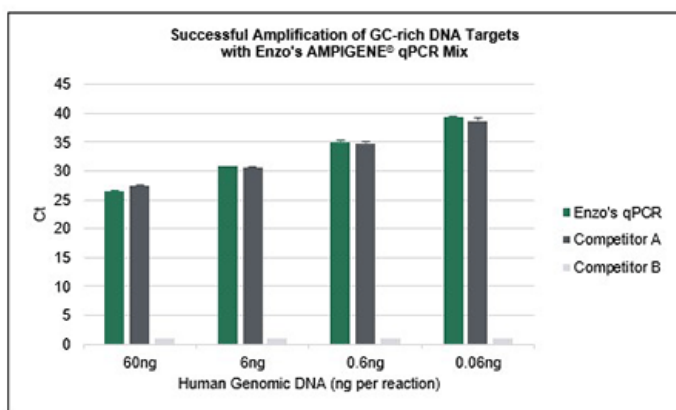


Figure 3. Amplification results for the GC-rich DNA target were obtained by targeting a human genomic region that contains 72.9% GC content. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system. The reactions with competitors' reagents were included and the performances were assessed.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

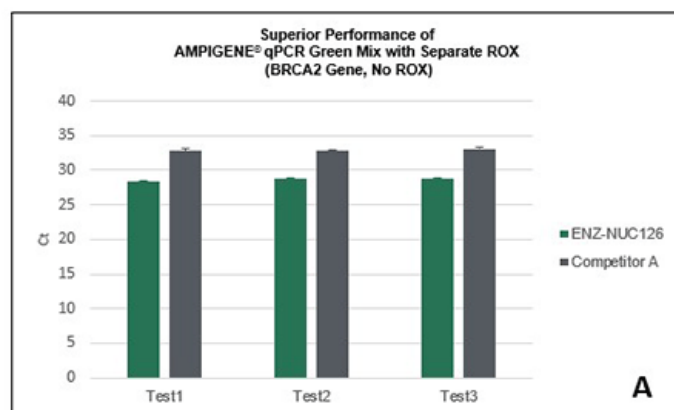


Figure 1A. Amplification results for the breast cancer gene 2 (BRCA2) were obtained from human genomic DNA samples in absence of ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product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

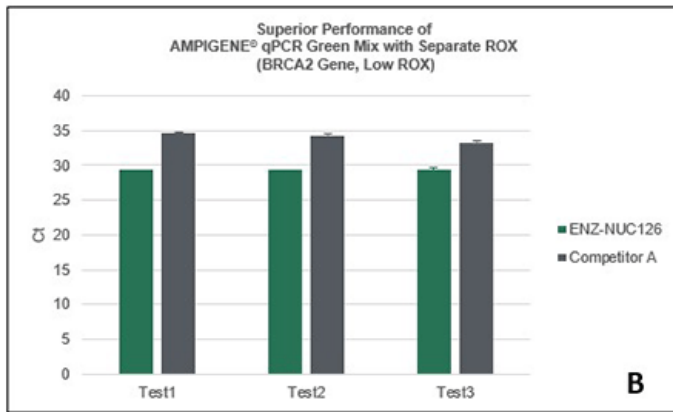


Figure 1B. Amplification results for the breast cancer gene 2 (BRCA2) were obtained from human genomic DNA samples with low ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were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

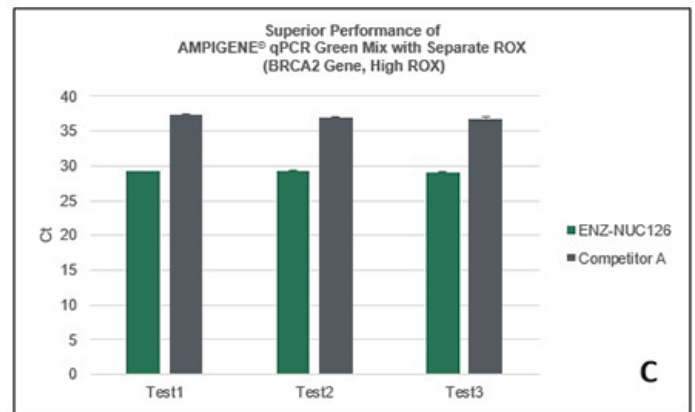


Figure 1C. Amplification results for the breast cancer gene 2 (BRCA2) were obtained from human genomic DNA samples with high ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were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

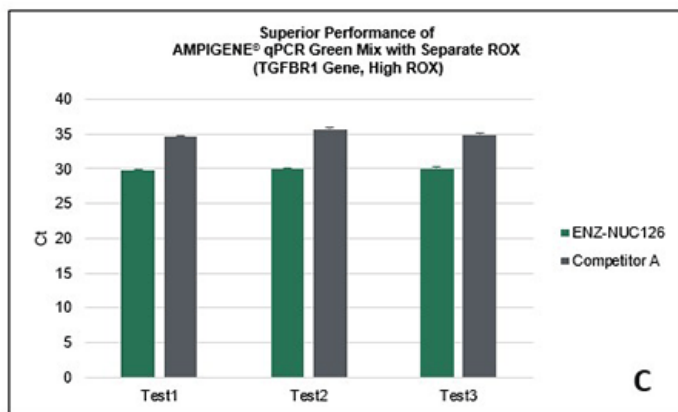


Figure 2C. Amplification results for the Transforming growth factor beta receptor 1 (TGFBR1) gene were obtained from human genomic DNA samples with high ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Green Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

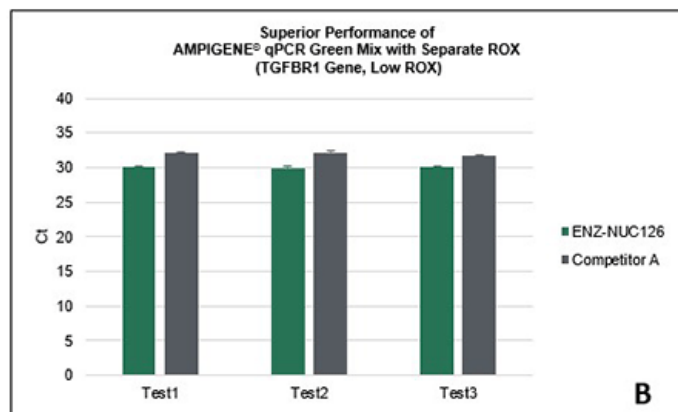


Figure 2B. Amplification results for the Transforming growth factor beta receptor 1 (TGFBR1) gene were obtained from human genomic DNA samples with low ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Green Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

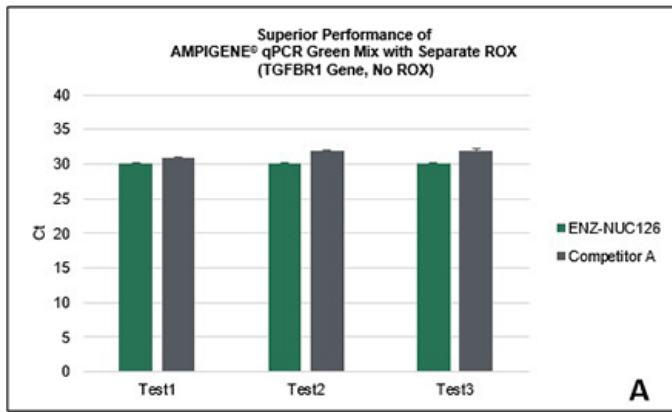


Figure 2A. Amplification results for the Transforming growth factor beta receptor 1 (TGFB1) gene were obtained from human genomic DNA samples in absence of ROX reference dye. Reactions were run in triplicate on a QuantStudio™ 5 real-time PCR system and each test was repeated three times. The reactions with competitor's product were included in each test and the performance was assessed. AMPIGENE® qPCR Green Mix with Separate ROX shows higher sensitivity in comparison with the competitor's product.

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application qPCR

Application Notes For dye-based qPCR

Contents 2X AMPIGENE[®] qPCR Green Mix (AMPIGENE[®] Hot Start
Taq DNA Polymerase, PCR green dye, 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)