

# CYGREEN<sup>®</sup> Nucleic Acid Dye

**Nucleic acid dye for gel staining and higher sensitivity qPCR**

CYGREEN<sup>®</sup> Nucleic Acid Dye is a DNA intercalating agent that is used to stain DNA. The DNA-dye complex emits a fluorescence spectra that makes it suitable for qPCR and gel staining applications.

CYGREEN<sup>®</sup> Nucleic Acid Dye has excitation/emission spectra similar to commonly used qPCR dyes and can be substituted in to increase sensitivity and reduce cost.

**Citations:** 10

[View Online »](#)

## Ordering Information

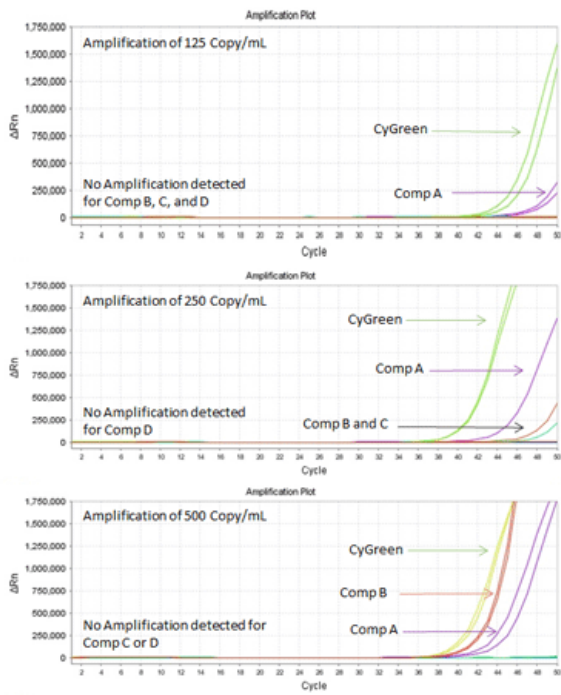
[Order Online »](#)

ENZ-GEN105-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

- Higher sensitivity allows detection of low target concentration
- Greater PCR efficiency for accurate and reproducible results
- Soluble in water for higher stability
- Compatible with various enzyme systems
- Suitable for commonly used rotor-type and plate-type qPCR platforms
- Validated for qPCR applications and gel staining
- Lower cost per assay compared to commonly used dyes



CYGREEN® Nucleic Acid Dye has greater sensitivity compared to competitor dyes

## Handling & Storage

Handling	Protect from light. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Reddish orange liquid.
Application	Gel staining, Melting curve, qPCR
Concentration	10,000x
Emission Maximum	515nm
Excitation Maximum	489nm
Formulation	Solution in DMSO.
Purity	≥93% (HPLC)



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)