



Glowing products for science

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

EvaGreen® Plus Dye, 20X in Water

EvaGreen® Plus Dye has an improved signal-to-noise compared to original EvaGreen® dye, for higher sensitivity in DNA amplification applications.



Product attributes

Excitation/Emission	487/525 nm (with DNA)
Concentration	20X

Product Description

EvaGreen® Plus Dye is an improved alternative to Biotium's original EvaGreen® Dye. EvaGreen® Plus Dye retains many of the essential benefits of the original patented EvaGreen® Dye, but has an improved signal-to-noise ratio that is more advantageous for a variety of DNA detection applications, including qPCR, digital PCR, and LAMP.

- Lower background & higher signal compared to original EvaGreen® Dye
- Better match for common qPCR instrument excitation sources
- Compatible with qPCR & HRM®
- Higher signal to noise offers advantages for digital PCR and isothermal applications
- Directly visualize amplification products in gel

Improvements over the Original

Similar to our original [EvaGreen® Dye](#), EvaGreen® Plus Dye is a green fluorescent nucleic acid dye that is essentially nonfluorescent by itself, but becomes highly fluorescent upon binding to dsDNA. EvaGreen® Plus Dye improves upon these properties with lower background fluorescence and increased brightness upon binding dsDNA. The dye has an excitation maximum that is a better fit for common qPCR instruments, resulting in improved sensitivity. As a result, EvaGreen® Plus Dye can give earlier Ct values and better signal discrimination. These properties offer potential advantages for digital PCR and isothermal amplification applications.

Why EvaGreen® Dye?

The unique properties of Biotium's EvaGreen® dyes make them particularly useful in quantitative real-time PCR applications. Relative to other green dyes, such as SYBR® Green I, both EvaGreen® dyes are generally less inhibitory toward PCR and less likely to cause nonspecific amplification. Both EvaGreen® dyes enable direct visualization of the PCR product on a gel using a UV transilluminator or blue light box without the need for another gel stain. Furthermore, EvaGreen® dyes can be used at a much higher dye concentration, resulting in more robust PCR signal.

EvaGreen® Plus Dye is provided at 20X in water. Also see our ready-to-use [Forget-Me-Not™ EvaGreen® qPCR Master Mix with ROX](#) as well as our [Forget-Me-Not™ EvaGreen® qPCR Master Mix with 2-Color Tracking](#) for sensitive, competitively-priced qPCR reagents. We also offer [Thiazole Green, 1000X in DMSO](#), which is identical to SYBR® Green I.

EvaGreen Dye and applications are covered by granted US and international patents. SYBR is a registered trademark of Thermo Fisher Scientific. HRM is a registered trademark of Idaho Technologies, Inc./BioFire Defense, LLC and its use may require a license.

References

Download list of curated [EvaGreen® Dye references](#).

This datasheet was generated on January 19, 2026 at 04:39:57 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/evagreen-plus-dye/>