

Helixyte Green™ Nucleic Acid Stain *200X* (Equivalent to PicoGreen®)

Ordering Information

Product Number: 17597 (1 mL in DMSO, 200X)
17598 (10 mL in DMSO, 200X)

Storage Conditions

Keep in -20°C. Avoid exposure to light

Biological Applications

Helixyte Green™ dsDNA stain is an ultra-sensitive fluorescent nucleic acid stain for quantitating double-stranded DNA (dsDNA) in solution. The Helixyte Green™ dsDNA stain has recently been used to quantitate PCR amplification yields in a method for direct cycle sequencing of PCR products. As little as 25 pg/mL of dsDNA (50 pg dsDNA in a 2 mL assay volume) with a standard spectra fluorimeter and 2.5 ng/mL dsDNA with a fluorescent microplate reader were detected with minimal effect in the presence of ssDNA, RNA, and free nucleotides. The assay is linear over three orders of magnitude and has little sequence dependence. It is ideal for accurately measuring DNA from many sources, including genomic DNA, viral DNA, miniprep DNA, or PCR.

Spectral Properties

Ex/Em = 502/523 nm

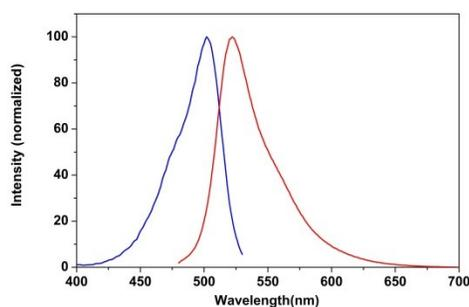


Figure 1. Excitation and emission spectra for the Helixyte Green™ dsDNA stain bound to DNA in PBS (pH 7.4).

Assay Protocol

The following protocol is an example for quantifying dsDNA with Helixyte Green™. Allow the Helixyte Green™ to warm to room temperature before opening the vial

Caution: No data are available addressing the mutagenicity or toxicity of Helixyte Green™ dsDNA stain. Because this reagent binds to nucleic acids, it should be treated as a potential mutagen and handled with appropriate care. The DMSO stock solution should be handled with particular caution as DMSO is known to facilitate the entry of organic molecules into tissues.

1. Preparing the Helixyte Green™ working solution:

- 1.1 Prepare an aqueous working solution of the Helixyte Green™ by making a 200-fold dilution of the concentrated DMSO solution in TE (10 mM Tris-HCl, 1 mM EDTA, pH 7.5–8.0). For example, add 50 µL Helixyte Green™ to 10 mL TE to prepare enough working solution to assay 100 samples in a 200 µL final volume. Protect the working solution from light by covering it with foil or placing it in the dark.

Note 1: We recommend preparing this solution in a plastic container rather than glass, as the dye may adsorb to glass surfaces.

Note 2: For best results, this solution should be used within a few hours of its preparation.

2. Prepare serial dilutions of dsDNA standard (0 to 3 ng/mL):

- 2.1 Prepare a 1 mg/mL stock solution of dsDNA (such as calf thymus DNA from Sigma) in ddH₂O.
- 2.2 Add 10 µL of 1 mg/mL dsDNA stock solution (from Step 2.1) to 998 µL TE buffer to have 10 µg/mL dsDNA solution, and then perform 1:10 and 1:2 serial dilutions to get 1000, 100, 50, 25, 12.5, 6.25, 3.125, and 0 ng/mL.
- 2.3 Add dsDNA standards and DNA containing test samples into a 96-well solid black microplate as described in Tables 1 and 2.

