

Product Information

NucBeacon™ Green Live nucleic acid stain

Catalog Number	Packaging Size
C013-1	1 mL

Storage upon receipt:

- -20°C
- Protect from light

Ex/Em: 500/530 nm, bound to DNA

Product Description

NucBeacon™ Green Live nucleic acid stain is cell-permeant nucleic acid stain that shows a large fluorescence enhancement upon binding nucleic acids. The NucBeacon™ Green Live nucleic acid stain can be used to stain RNA and DNA in both live and dead eukaryotic cells, as well as in Gram-positive and Gram-negative bacteria. The novel NucBeacon™ Green Live nucleic acid stain has several important characteristics:

- Permeability to virtually all cell membranes, including mammalian cells and bacteria.
- High molar absorptivity, with extinction coefficients $>60,000 \text{ cm}^{-1} \text{ M}^{-1}$ at visible absorption maxima.
- Extremely low intrinsic fluorescence, with quantum yields typically <0.01 when not bound to nucleic acids.
- Quantum yields that are typically >0.4 when bound to nucleic acids.

The NucBeacon™ Green Live nucleic acid stain is compatible with a variety of fluorescence-based instruments that use laser excitation or a conventional broadband illumination source (e.g., mercury- and xenon-arc lamps). The NucBeacon™ Green Live nucleic acid stain has been used in diverse applications from staining DNA spotted on microarrays to staining live and fixed cells.

Experimental Guidelines

We suggest broad ranges of staining concentrations, based on our laboratory experience or published methods, to provide a starting point for experiments. These conditions require adjustment for each cell type and experimental system.

Use plastic tubes when diluting the NucBeacon™ Green Live nucleic acid stain, because the diluted stain adheres to glass.

In general, the best results are obtained in buffers that do not contain phosphate. When preparing other solutions, note that residual detergent on plastic or glassware may also affect real or apparent staining of many cells or organisms, causing brightly stained material to appear in solutions with or without cells present. Wash all labware in mild detergent and rinse

with hot tap water followed by several rinses with deionized water.

Adherent cells in culture may be stained *in situ* on coverslips. Pellet cells in suspension by centrifugation and resuspend in buffered salt solution or water. Add the NucBeacon™ Green Live nucleic acid stain using the concentrations listed in Table 1 as a guide. In initial experiments, it may be best to try several dye concentrations over the entire suggested range to determine the concentration that yields optimal staining. Be aware that growth medium, cell density, the presence of other cell types, and other factors may influence staining. Stained eukaryotic cells generally show diffuse cytoplasmic staining as well as nuclear staining. Particularly intense staining of intranuclear bodies is frequently observed. Because the NucBeacon™ Green Live nucleic acid stain is cell permeant and contain a net positive charge at neutral pH, they may also stain mitochondria. Staining of live yeast is primarily mitochondrial.

The NucBeacon™ Green Live nucleic acid stain has proven to be useful for staining DNA on microarrays for quality control purposes.

Table 1. Recommended conditions for staining cells with the NucBeacon™ Green Live nucleic acid stain.

Application	Dye Concentration	Staining Conditions
Bacteria	50 nM–20 μM	Incubate for 1–30 minutes.
Eukaryotic cells	10 nM–5 μM	Incubate for 10–120 minutes
Microarrays	50 nM in TE buffer	Incubate for 5 minutes, rinse and then dry.