

ab138904

Nuclear Green LCS1

Instructions for Use

The labelling of DNA in cells

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Biological Applications	3
2. Spectral Properties	3
3. Sample Protocol for Cell Staining	4

1. Biological Applications

ab138904 Nuclear Green LCS1 is a non-fluorescent, DNA-selective and cell-permeant dye for analyzing DNA content in living cells. The Nuclear Green LCS1 has its green fluorescence significantly enhanced upon binding to double-stranded DNA. It can be used in fluorescence imaging, microplate and flow cytometry applications. This DNA-binding dye might be used for multicolor analysis of live cells.

2. Spectral Properties

Ex/Em = 503/526 nm

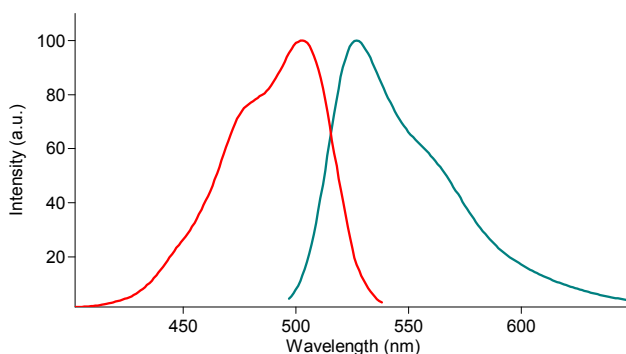


Figure 1. Excitation and emission spectra for the Nuclear Green LCS1 bound to DNA in PBS (pH 7.4).

3. Sample Protocol for Cell Staining

Caution: The following protocol can be adapted for most cell types. Growth medium, cell density, the presence of other cell types and factors may influence staining. Residual detergent on glassware may also affect staining of many organisms, and cause brightly stained material to appear in solutions with or without cells present.

1. Add Nuclear Green LCS1 (2 to 10 μM) into the cells (either suspension or adherent cells), and stain the cells for 15 to 60 minutes. In initial experiments, it may be best to try several dye concentrations to determine the optimal concentration that yields the desired result. High dye concentration tends to cause nonspecific staining of other cellular structures.
2. Directly analyze the cellular staining with a fluorescence microscope, a fluorescence microplate reader, or flow cytometer.

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “contact us” on www.abcam.com for the phone number for your region).

UK, EU and ROW

Email: technical@abcam.com

Tel: +44 (0)1223 696000

www.abcam.com

US, Canada and Latin America

Email: us.technical@abcam.com

Tel: 888-77-ABCAM (22226)

www.abcam.com

China and Asia Pacific

Email: hk.technical@abcam.com

Tel: 108008523689 (中國聯通)

www.abcam.cn

Japan

Email: technical@abcam.co.jp

Tel: +81-(0)3-6231-0940

www.abcam.co.jp