

## Nuclear Red™ LCS2

### Ordering Information

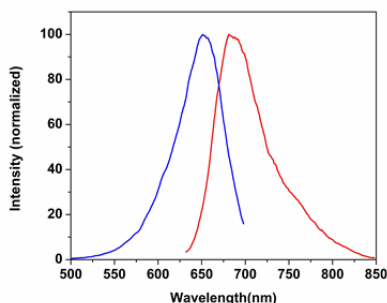
Product Number: 17545 (5 mM)

### Storage Conditions

Keep at -20 °C and avoid exposure to light

### Spectral Properties

Ex/Em = 651/681 nm



**Figure 1.** Excitation and emission spectra for the Nuclear Red™ LCS2 bound to DNA in PBS (pH 7.4).

### Biological Applications

Our Nuclear Red™ LCS2 is a fluorogenic, DNA-selective and cell-permeant dye for analyzing DNA content in living cells. The Nuclear Red™ LCS2 has its red fluorescence significantly enhanced upon binding to 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 with the Cy5 filter sets.

### 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 2 to 10  $\mu$ M 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.

**Disclaimer:** This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.