

## Product information: SPY595-DNA (SC301)

Live Cell Fluorogenic DNA Labelling Probe

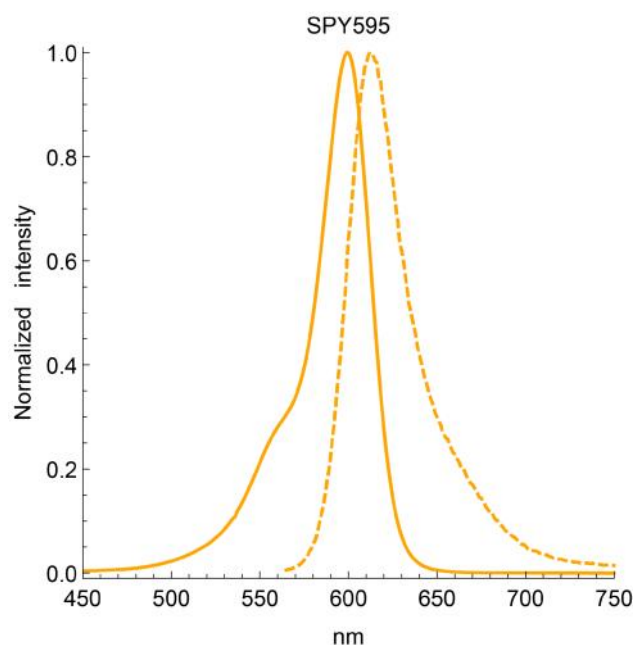
### Introduction

SPY595-DNA is a bright red & non toxic live cell nuclear stain based on our SPY™ dyes series. Its optimized structure allows quick labeling of DNA in live & fixed cells with high specificity and very low background. SPY595-DNA stains the nuclei of live or fixed cells without the need for genetic manipulation or overexpression of fluorescent proteins. Its emission in the red minimizes phototoxicity and sample autofluorescence. SPY595-DNA enables multicolor imaging with SPY505, SPY555, SPY650, SPY700, SiR and GFP. SPY595-DNA can be imaged with standard Alexa594™ or Texas Red filtersets. It can be used for widefield, confocal, SIM or STED imaging in living or fixed cells and tissue. Contains 1 vial of SPY595-DNA (lyophilized).



### Probe Properties

<b>Absorbance maximum</b> <sub>abs</sub>	599 nm
<b>Fluorescence maximum</b> <sub>fl</sub>	615 nm
<b>Works on fixed cells?</b>	yes, PFA and methanol
<b>Probe quantity</b>	100 stainings*
<b>Fluorescence lifetime</b>	n.d.
<b>STED depletion wavelength</b>	775 nm
<b>Shipping</b>	room temperature
<b>Storage</b>	-20°C



### Storage & Handling

Store the probe at -20°C or below upon receipt. The lyophilized probe is stable for >1 week at room temperature and for >12 months at -20°C. Reconstitute SPY595-DNA using anhydrous DMSO. We recommend to use newly or freshly opened and anhydrous DMSO to prepare the 1000x stock solution. In contact with air and moisture, DMSO produces decay products which can strongly reduce the shelf life of the probe in solution, even at -20°C. Keep the 1000x stock solution of the compound below -20°C after use. Vials should be allowed to warm to room temperature before opening. When reconstituted and stored properly, the 1000x stock solution is stable for 3 months. Note: DMSO solutions should be handled with particular caution as DMSO is known to facilitate the entry of organic molecules into tissues. Dispose of these reagents in compliance with all pertaining local regulations.

### Labelling Protocol

**Note:** This protocol was optimized using HeLa cells adhering to coverslips and has been confirmed in other common cell lines. Recommendations in this protocol should be used as a starting point, and optimal labeling conditions for each cell type should be determined empirically. SPY595-DNA is based on the DNA minor groove binding molecule bisbenzimidazole. It may modify DNA metabolism in living cells at high doses. Therefore the recommended staining dilution is 1000 fold or more if long term (>12h) imaging experiments are planned. For all other purposes, 1000 fold dilution SPY595-DNA for staining is recommended.

**1. Prepare 1000x stock solution.** Add 50  $\mu$ L of anhydrous DMSO to the SPY595-DNA vial to prepare the 1000x stock solution. We recommend to use newly or freshly opened and anhydrous DMSO to prepare the 1000x stock solution. In contact with air and moisture, DMSO produces decay products which can strongly reduce the shelf life of the probe in solution, even at  $-20^{\circ}\text{C}$ . At this stage, the solution can be colored or not, this has no influence on the performance of the probe. After use, this solution should be stored at  $-20^{\circ}\text{C}$  or below. Do not divide the 1000x stock solution into small aliquots, they will decay faster and the probe is not altered by multiple freeze-thaw cycles. When stored properly, this stock solution is stable for 3 months.

**2. Prepare the staining solution.** Dilute SPY595-DNA to 1x in your usual cell culture medium (e.g. DMEM + 10% fetal bovine serum) and vortex briefly. If the dilution is not performed in a single step, please use DMSO to prepare the intermediate dilution as using aqueous buffers to prepare the intermediate dilution will lead to the formation of probe aggregates. Proceed quickly to step 3. Since staining efficiency can depend on the cell line, it is recommended to stain cells with 1000x dilution at the first attempt and then optimize the SPY595-DNA dilution factor in further experiments until an optimal staining is achieved (see labelling concentration & incubation time table below). Use only freshly made staining solution, and do not use it multiple times.

**3. Cell preparation and staining.** Grow cells on coverslips, glass bottom dish or glass bottom multi-well plates as usual. When cells have reached the desired density, replace the culture medium by the **staining solution** freshly prepared under step 2 ensuring that all the cells are covered with the solution. Place the cells in the incubator at  $37^{\circ}\text{C}$  in a humidified atmosphere containing 5%  $\text{CO}_2$  and observe the following table to determine labelling time as a function of probe concentration:

Dilution factor	suggested labelling time (h)**
1000 or less	1
2000	2
>2000	4

**Note: SPY595-DNA stains DNA in paraformaldehyde (PFA) and methanol fixed cells.**

**4. Cell imaging.** Imaging of SPY595-DNA is best performed using standard Alexa594<sup>TM</sup> or Texas Red settings. After labelling, the live cells can be immediately imaged without the need for washing steps. Optionally, a simple washing step consisting of replacing once the labelling solution by fresh culture medium which does not contain the probe may improve the signal to noise ratio. If time lapse imaging is performed, it is recommended to keep the probe in the imaging medium during the whole experiment to get a constant signal. If the cells were washed before imaging, the staining will last for a few hours. Please note that SPY595-DNA may be excited by 360-390 nm light and produce some fluorescence at ca. 450 nm due to fluorescence of the DNA binding moiety of the probe.

\* Based on the following conditions: 0.5 ml staining solution / staining experiment with 1x probe concentration. The number of staining experiments can be further increased by reducing volume or probe concentration.

\*\* These labelling times were determined for HeLa cells and may differ depending on the cell line used.

Spirochrome products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply. Spirochrome products and product applications are covered by patents and patents pending. SPY is a registered trademark

**Limited Use Label License:** For research use only. Not intended for any animal or human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Spirochrome will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Spirochrome is willing to accept return of the unused product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Spirochrome: Spirochrome AG, Postfach 213, 8620 Stein am Rhein, Switzerland, Email: [info@spirochrome.com](mailto:info@spirochrome.com)