

## Product information: SiR700-tubulin Kit (CY-SC014)

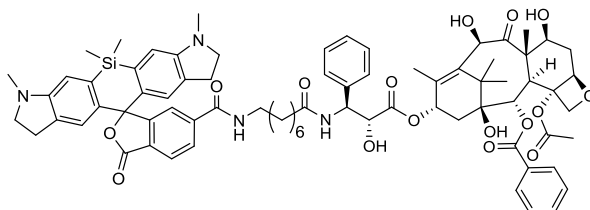
Live Cell Fluorogenic Microtubule Labelling Probe

### Introduction

SiR700-tubulin is based on the silicon rhodamine (SiR) fluorophore analogue SiR700 and the microtubule binding drug Docetaxel. SiR700-tubulin allows the labelling of microtubules in live cells with high specificity and low background<sup>1</sup>. The key features of SiR700-tubulin are i) far-red absorption and emission wavelengths, ii) cell permeability, iii) fluorogenic character and iv) compatibility with superresolution microscopy (STED & SIM). The unprecedented combination of those properties in a single probe put SiR700-tubulin at the leading edge of excellence.

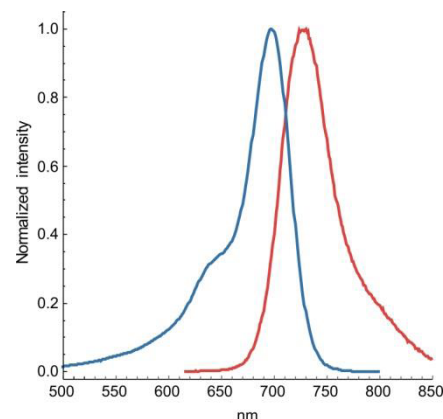
### Physical properties

abs	689 nm
Em	716 nm
$\epsilon_{689 \text{ nm}}$	$1.0 \cdot 10^5 \text{ mol}^{-1} \cdot \text{cm}^{-1}$
MW	1327.6 g/mol
MF	$\text{C}_{75}\text{H}_{86}\text{N}_4\text{O}_{16}\text{Si}$



### Storage & Handling

Store the compound below  $-20^\circ\text{C}$  upon receipt. Prepare solutions of the compound using anhydrous DMSO. Keep solutions of the compound below  $-20^\circ\text{C}$  after use. Vials should be allowed to warm to room temperature before opening. When stored properly, the compound should be stable for several 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.



**Kit contents:** 35 nmol SiR700-tubulin and 1  $\mu\text{mol}$  verapamil

### Labelling Protocol

**Note:** This protocol was optimized using human fibroblast cells adhering to coverslips and has been confirmed in other common cell lines. Recommendations for experimental protocols should be used as a starting point, and optimal labeling conditions for each cell type should be determined empirically. SiR700-actin is based on the actin stabilizing drug jasplakinolide. It can therefore modify tubulin dynamics in living cells. Whereas interphase cells were only little affected by the probe, concentrations above 100 nM of SiR700-tubulin led to mitotic duration increase in cultured HeLa cells<sup>1</sup>. If long term imaging experiments are planned where tubulin dynamics are critical, we recommend to keep the concentration of SiR700-tubulin equal or below 100 nM. For other purposes, using 0.5-1  $\mu\text{M}$  SiR700-tubulin for staining is recommended.

**Prepare 1 mM stock solution.** Dissolve the content of the vial of SiR700-tubulin in 35  $\mu\text{L}$  of anhydrous DMSO to make a 1 mM stock solution. This solution should be stored at  $-20^\circ\text{C}$  or below. **Do not divide the solution into small aliquots**, they will decay faster and the compound is not altered by multiple freeze-thaw cycles. When stored properly, this stock solution should be stable for three months or more. If the concentration of the stock solution needs to be accurately determined, dilute 1  $\mu\text{L}$  of 1 mM stock solution in 99  $\mu\text{L}$  of PBS containing 0.2 % SDS. After 15 minutes at room temperature, measure the absorbance at 652 nm. Calculate the concentration using the extinction coefficient given above.

**Prepare staining solution.** Dilute SiR700-tubulin to the desired concentration in cell culture medium (e.g. DMEM + 10% fetal bovine serum) and vortex briefly. Since staining efficiency can depend on the cell line, it is recommended to stain cells with 1  $\mu\text{M}$  at the first attempt to quickly obtain a strong staining and then reduce the SiR700-tubulin concentration in further experiments until an optimal staining is achieved (see labelling concentration & incubation time table below). Some cell lines might express high levels of efflux pumps and may therefore be poorly stained by SiR700-tubulin. The addition of 10  $\mu\text{M}$  verapamil, a broad spectrum efflux pump inhibitor, in the staining solution usually greatly improves the staining. Use only freshly made staining solution and do not use it multiple times.

**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** ensuring that all the cells are covered with solution. Place the cells in the incubator at 37°C in a humidified atmosphere containing 5% CO<sub>2</sub> and observe the following table to determine labelling time as a function of probe concentration:

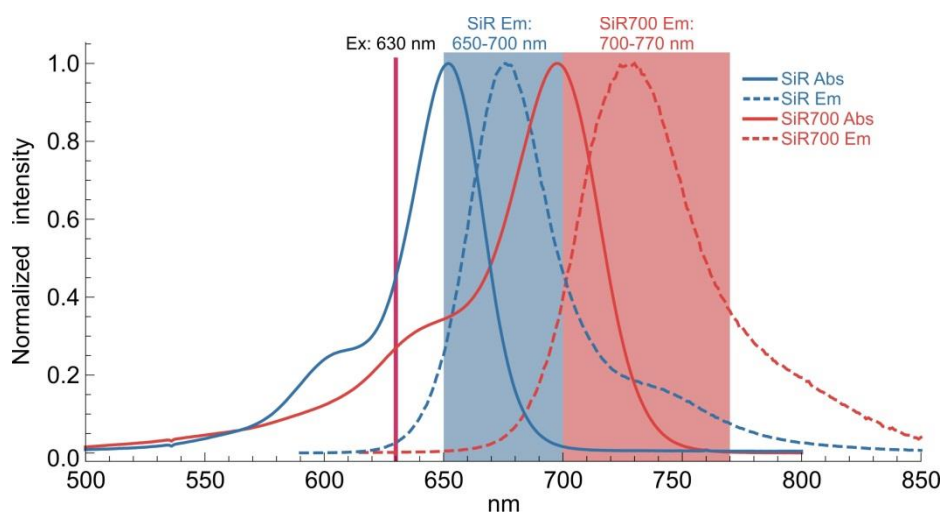
probe concentration (nM)	suggested labelling time (h)*
> 1000	0.5 - 1
500	3 - 4
200	4 - 6
< 100	6 - 12

\* these labelling times were determined for human fibroblasts and may differ depending on the cell line used.

**Note: SiR700-tubulin does not stain paraformaldehyde (PFA) and methanol fixed cells as these fixation methods alter the probe binding site on microtubules.**

**Cell imaging.** Imaging of SiR700-tubulin can be performed using standard Cy5 settings, however recording fluorescence emission between 700 and 800 nm provides a better signal. After labelling, the live cells can be immediately imaged without the need for washing steps. Optionally, a simple washing step consisting of replacing 1-3 times the labelling solution by fresh culture medium which does not contain the probe usually improves the signal to noise ratio. If time lapse imaging is performed, it is recommended to keep the concentration of probe equal or below 100 nM during the whole experiment to get a constant signal and to avoid interference of the probe with tubulin dynamics (reduced cell proliferation). If cells were washed before imaging, the staining will last for a few hours.

**Dual colour imaging.** SiR700-tubulin can be used together with spirochrome SiR-probes to perform dual colour imaging. SiR700-tubulin and the second SiR-probe (e.g. SiR-actin, SiR-DNA or SiR-lysosome) can be added simultaneously to the cells using the standard protocol above. A washing step is recommended to obtain the best signal to background ratio. Imaging is best performed using the following parameters: excitation using 630 to 640 nm light for both SiR700 and SiR; emission 650-700 nm (SiR channel) and 700-770 nm (SiR700 channel). A small bleed through of SiR into the SiR700 channel and vice versa may be observed. For more information and examples of dual colour imaging please visit our website [www.spirochrome.com/dualcolour](http://www.spirochrome.com/dualcolour).



## References:

1. Fluorogenic probes for live-cell imaging of the cytoskeleton, G. Lukinavičius et al., *Nature Methods*, 11, 731–733 (2014)
2. Fluorogenic probes for multicolor imaging in living cells, G. Lukinavičius et al., *JACS*, (2016)

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.

**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)