

**Tubulite™ Red \*Cell-Permeable\***

 Catalog number: 23718  
 Unit size: 100 Slides

Component	Storage	Amount (Cat No. 23718)
Tubulite™ Red	Freeze (< -15 °C), Minimize light exposure	100 Slides

**OVERVIEW**

Tubulite™ Red is a robust tool for fluorescent visualization of tubulin structures in live cells. This cell-permeable probe enables real-time imaging of tubulin dynamics, bypassing the need for cell fixation and allowing continuous assessment of tubulin polymerization and organization. The probe's red fluorescence and high cellular permeability facilitate its use in multiplexed imaging applications, including co-staining with GFP and nuclear dyes such as DAPI. Upon cellular entry, Tubulite™ Red is hydrolyzed by intracellular esterases, removing its lipophilic blocking group to produce a charged, membrane-impermeant form that is well-retained within cells. Tubulite™ Red is not generally recommended for studies of dynamic cellular processes, as it may interfere with key cellular functions, including tubulin activity and cell division. Additionally, it is unsuitable for fixed-cell imaging, as fluorescence is lost during cell fixation.

**AT A GLANCE**
**Protocol Summary**

1. Prepare cells with test compounds at a density of  $5 \times 10^5$  to  $1 \times 10^6$  cells/mL
2. Prepare and add Tubulite™ Red working solution to cells
3. Incubate at 37°C for 60 minutes.
4. Use a fluorescence microscope with a Cy3 filter set to measure fluorescence intensity.

**Note:** This protocol is intended as a general guide and should be adjusted to meet your specific requirements.

**KEY PARAMETERS**
**Fluorescence microscope**

Emission	Cy3 filter set
Excitation	Cy3 filter set
Recommended plate	Black wall/clear bottom

**CELL PREPARATION**
**Prepare Cells**

1. For each sample, prepare cells in 1 mL of warm medium or buffer at a density of  $5 \times 10^5$  to  $1 \times 10^6$  cells/mL.

**Note:** Each cell line should be individually assessed to determine its optimal cell density.

**PREPARATION OF STOCK SOLUTIONS**

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles

**Tubulite™ Red Stock Solution (100X)**

1. Add 100 µL of DMSO (not provided) to the Tubulite™ Red vial and mix thoroughly.

**Note:** Aliquot and store any unused Tubulite™ Red stock solution at -20°C. Avoid multiple freeze-thaw cycles.

**PREPARATION OF WORKING SOLUTION**
**Tubulite™ Red Working Solution (1X)**

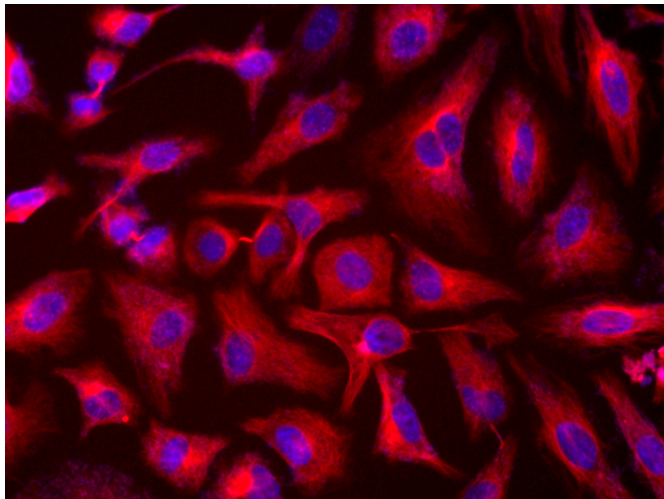
1. To prepare a 1X Tubulite™ Red working solution, add 10 µL of Tubulite™ Red stock solution and 1 mM ReadiUse™ Probenecid (AAT Cat# 20061, not provided) to 1 mL of HHBS [Hanks' Buffer with 20 mM Hepes] (AAT Cat# 20011, not provided) or your preferred buffer. Mix thoroughly.

**Note:** For best results, we recommend preparing a fresh Tubulite™ Red working solution each time you use it. The working solution remains stable for a few hours.

**Note:** The calculations above assume that 100 µL of working solution will be used for each slide or sample. Please adjust the concentrations of Tubulite™ Red and ReadiUse™ probenecid to match your specific staining and cell conditions.

**SAMPLE EXPERIMENTAL PROTOCOL**

1. Prepare cell samples as needed.
  2. Remove the cell growth medium, then wash the cells with PBS (not included) or another buffer of your choice.
  3. Add 100 µL of Tubulite™ Red working solution to the cells, then place them in a 37°C incubator for 60 minutes.
- Note:** The incubation time will vary based on the specific cell type and cell concentration. Be sure to adjust the incubation time to best suit each experiment.
4. Remove the working solution and wash the cells twice with PBS or a preferred buffer.
  5. Cover the cells with HHBS (or a buffer of your choice) containing 1 mM probenecid. Then, use a fluorescence microscope equipped with a Cy3 filter set to monitor the fluorescence intensity.



**Figure 1.** Detection of Tubulin in HeLa Cells: HeLa cells were co-labeled with Tublite™ Red and DAPI (AAT Bioquest Cat# 17507) for 60 minutes in a 37°C incubator with 5% CO<sub>2</sub>. Fluorescence intensity was measured using a fluorescence microscope with the Cy3 channel.

**DISCLAIMER**

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.