

ReadiCleave™ AML Cleavage Buffer *5X Aqueous Solution*

Catalog number: 7510

Unit size: 1 mL

Component	Storage	Amount (Cat No. 7510)
ReadiCleave™ AML Cleavage Buffer *5X Aqueous Solution*	Freeze (< -15 °C), Minimize light exposure	1 mL

OVERVIEW

ReadiCleave™ AML Cleavage Buffer provides an optimized formula to effectively remove the fluorescence tags from our ReadiCleave™ AML-labeled biological targets when needed. The removal of the fluorescent ReadiCleave™ AML tags may make the fluorescence imaging more compatible with the downstream analysis of a biological target by other techniques, e.g., MS analysis. Using ReadiCleave™ AML Cleavage Buffer alongside ReadiCleave™ Styramides enables multiplexed fluorescence imaging applications in cells and tissues. ReadiCleave™ Styramides are the new iteration of our PSA products that add the reversible capability to chemically remove the PSA staining on tissue or in cells (if desired) while preserving the integrity of tissue samples.

PREPARATION OF WORKING SOLUTION
ReadiCleave™ AML Cleavage Buffer Working Solution (1X)

- To prepare a 1X working solution, add 200 µL of ReadiCleave™ AML Cleavage Buffer into 800 µL of ddH₂O, and mix thoroughly.

Note: For optimal results, use this solution within a few hours of preparation.

SAMPLE EXPERIMENTAL PROTOCOL

- Remove the coverslip by briefly dipping the slides into PBS. Then, wash the tissue sample with PBS to eliminate any remaining mounting medium.
- Add 100 µL of ReadiCleave™ AML Cleavage Buffer working solution to the tissue or cell samples.

Note: Add a sufficient amount of ReadiCleave™ AML Cleavage Buffer working solution to ensure that the samples are fully submerged.
- Heat the samples at 60°C for 60 minutes.
- Remove the ReadiCleave™ AML Cleavage Buffer working solution and briefly rinse the samples with PBST.
- Reprocess the tissue samples beginning with the Antigen Retrieval step in your IHC staining protocol.

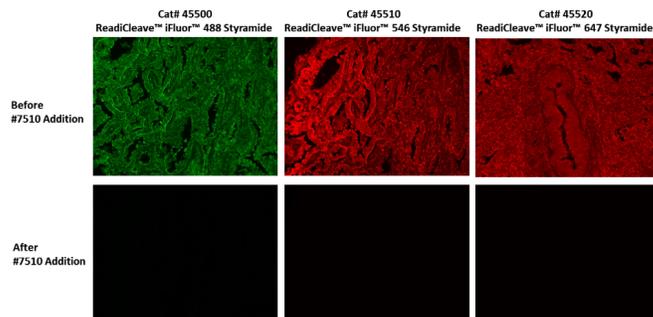
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. Fluorescence imaging for the detection of EpCAM in lung adenocarcinoma was conducted using ReadiCleave™ iFluor® 488 Styramide (Cat# 45500), ReadiCleave™ iFluor® 546 Styramide (Cat# 45510), and ReadiCleave™ iFluor® 647 Styramide (Cat# 45520). Images were captured both before and after treatment with ReadiCleave™ AML Cleavage Buffer, providing comparative insights into the effectiveness of the cleavage buffer treatment.

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 if you have any questions.