

## Tetracysteine REASH Reagent \*2 mM DMSO Solution\*

Catalog number: 22332  
Unit size: 100 Tests

| Component     | Storage                                    | Amount (Cat No. 22332) |
|---------------|--------------------------------------------|------------------------|
| REASH Reagent | Freeze (< -15 °C), Minimize light exposure | 100 Tests              |

### OVERVIEW

Tetracysteine REASH Reagent is a resorufin-derived biarsenical probe modified with two arsenic atoms positioned to interact with tetracysteine-containing protein tags. The biarsenical labeling technology is based on the high-affinity interaction between arsenic groups and thiols. Tetracysteine REASH Reagent binds recombinant proteins engineered with a tetracysteine sequence, typically a Cys-Cys-X1-X2-Cys-Cys motif. When REASH binds to the tetracysteine sequence, its biarsenical group reacts rapidly with the Cys-Cys moiety, generating a strong red fluorescent signal. This compact tag-based labeling approach is useful for studying protein localization, turnover, trafficking, receptor signaling, and internalization.

### AT A GLANCE

#### Protocol summary

1. Prepare cells
2. Prepare REASH Reagent working solution
3. Incubate the cells with REASH Reagent working solution for 15-60 minutes
4. Image the cells using the Cy3 filter set

### KEY PARAMETERS

#### Fluorescence microscope

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

### PREPARATION OF WORKING SOLUTION

#### REASH Reagent working solution (1X)

Dilute the 2mM stock solution at 1:800 in an appropriate buffer such as serum-free or low-level serum (~1%) medium, HHBS, or Opti-MEM® medium to make a 2.5 µM (1X) labeling solution and mix well.

**Note:** Make the working solution just before use.

**Note:** For cells transduced with lentivirus, a 0.5X working solution may be optimal. Depending on the levels of specific and background fluorescent signal, you can optimize the working solution to better visualize your labeled protein. We recommend trying a concentration range of 0.4 to 4X working solution.

**Table 1.** The following table is the suggested volume of labeling solution to use for different tissue culture formats.

| Plate  | 96-well | 48-well | 24-well | 12-well | 6-well |
|--------|---------|---------|---------|---------|--------|
| Volume | 100 µL  | 150 µL  | 250 µL  | 500 µL  | 1 mL   |

### SAMPLE EXPERIMENTAL PROTOCOL

1. Prepare cells as desired.
2. Remove the growth medium from the cells and wash cells once

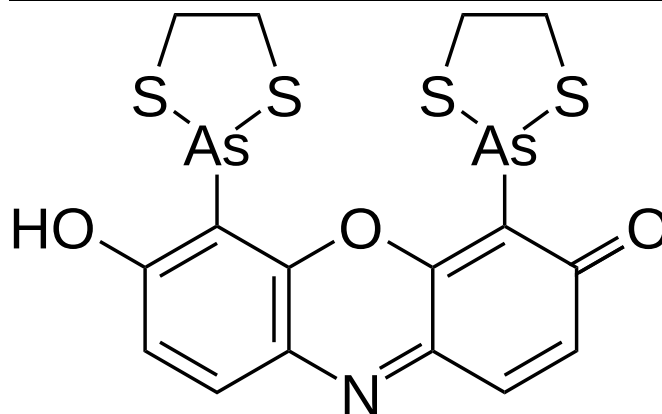
with an appropriate medium.

3. Add the appropriate amount of 1X REASH Reagent working solution to each well (See table for the appropriate volume).

**Note:** Discard any unused 1X working solution. Do not reuse the 1X working solution.

4. Incubate the cells at room temperature for 15-60 minutes, protected from light.
5. Wash the cells with 250 µM BAL in serum free medium or buffer of your choice 2 times.
6. Image your labeled protein using a fluorescence microscope with a Cy3 filter set.

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Chemical structure for REASH Reagent.

### 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.