

Thiolite™ Green

Ordering Information	Storage Conditions
Product Number: 21508 (5 mg)	Store at -20 °C, desiccated and protected from light Expiration date is 12 months from the date of receipt

Introduction

Thiolite™ Green is one of the most sensitive sensors for measuring thiol compounds. It gives a green fluorescent adduct upon reacting with thiol compounds (such as cysteine). It can be used to quantify the number of cysteines on a protein. When we use it to measure glutathione fluorimetrically, it has >200 fold fluorescence enhancement upon reaction with thiol-containing compounds.

Chemical and Physical Properties

Molecular Weight: ~400

Solvent: dimethylsulfoxide (DMSO)

Spectral Properties: Excitation = 510 nm; Emission = 524 nm

Assay Protocol with Thiolite™ Green in a 96-well Plate

Brief Summary

Prepare Thiolite™ Green working solution (50 µL) → Add GSH standards or test samples (50 µL)
→ Incubate at room temperature for 10 min-1 hr → Read fluorescence intensity at Ex/Em = 490/525 nm

Note: Following is our recommended protocol for thiol assay in solution. This protocol only provides a guideline, and should be modified according to your specific needs.

1. Prepare Thiolite Green™ working solution:

- 1.1 Prepare a 10 to 25 mM stock solution of Thiolite™ Green in high-quality, anhydrous DMSO. The stock solution should be used promptly; any remaining solution need be aliquoted and frozen at -20 °C.

Note: The unused Thiolite Green stock solution should be divided into single use aliquots and stored at -20°C, protected from light.

- 1.2 Prepare a 2X Thiolite™ Green working solution: On the day of the experiment, either dissolve Thiolite™ Green in DMSO or thaw an aliquot of the Thiolite™ Green stock solution to room temperature. Prepare a 2X working solution at the final concentration ranging from 100 to 250 µM in 20 mM Hepes buffer or buffer of your choice, pH 7. It is recommended to use Thiolite™ Green at the final concentration ranging from 50 to 100 µM to measure Thiol concentration in solution.

2. Run GSH Assay in supernatants:

- 2.1 Add 50 µL of 2X Thiolite™ Green working solution (from Step 1.2) to each well of the GSH standard, blank control, and test samples to make the total GSH assay volume of 100 µL/well.

Note: For a 384-well plate, add 25 µL of sample and 25 µL of GSH reaction mixture into each well.

- 2.2 Incubate the reaction at room temperature for 10 to 60 minutes, protected from light.

- 2.3 Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 490/525 nm.

2.4 The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with the thiol reaction.

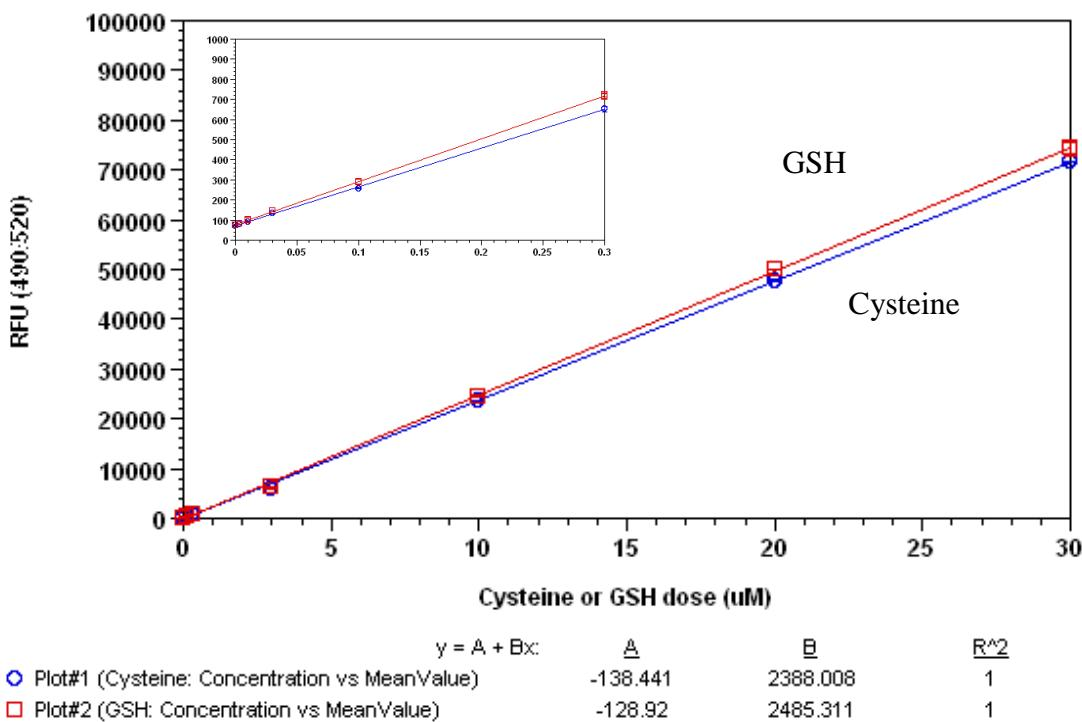


Figure 1. GSH and cysteine dose response was measured with Thiolite™ Green on a 96-well black plate using a NOVOStar microplate reader (BMG Labtech). As low as 10 nM (1 pmol/well) of GSH or cysteine can be detected with 10 minutes incubation time (n=3). The insert shows the low levels of thiol detection.

Disclaimer: This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.