

# Ellman's Reagent

22582

0311.3

Number	Description
22582	<b>Ellman's Reagent</b> , 5,5'-Dithio- <i>bis</i> -(2-nitrobenzoic acid), 5g Molecular Weight: 396.35 CAS #: 69-78-3

**Storage:** Store product protected from moisture at 4°C. Product is shipped at ambient temperature.

## Table of Contents

Introduction .....	1
Procedure for Quantitating Sulfhydryl Groups Using a Cysteine Standard.....	2
Procedure for Quantitating Sulfhydryl Groups Based on Molar Absorptivity.....	2
Troubleshooting.....	3
Related Thermo Scientific Products .....	3
Additional Information .....	4
Cited References.....	4

## Introduction

In 1959 Ellman<sup>1</sup> introduced 5,5'-dithio-*bis*-(2-nitrobenzoic acid), also known as DTNB, as a versatile water-soluble compound for quantitating free sulfhydryl groups in solution (Figure 1, see Additional Information section). A solution of this compound produces a measurable yellow-colored product when it reacts with sulfhydryls. Consequently, Thermo Scientific Ellman's Reagent is very useful as a sulfhydryl assay reagent because of its specificity for -SH groups at neutral pH, high molar extinction coefficient and short reaction time.

DTNB reacts with a free sulfhydryl group to yield a mixed disulfide and 2-nitro-5-thiobenzoic acid (TNB; see Figure 2). The target of DTNB in this reaction is the conjugate base ( $R-S^-$ ) of a free sulfhydryl group. Therefore, the rate of this reaction is dependent on several factors:<sup>2</sup> 1) the reaction pH, 2) the pKa' of the sulfhydryl and 3) steric and electrostatic effects. TNB is the "colored" species produced in this reaction and has a high molar extinction coefficient in the visible range. The molar extinction coefficient of TNB was originally reported to be  $13,600M^{-1}cm^{-1}$  at 412nm and pH 8.0.<sup>1</sup> Consequently, this value has been cited frequently in the literature. Later studies have shown, however, that the molar extinction coefficient is more accurately reflected by a value of  $14,150M^{-1}cm^{-1}$  at 412nm.<sup>2,3</sup> The extinction of TNB is not affected by changes in pH between 7.6 and 8.6.<sup>3</sup> However, the extinction of TNB is different in other solvents (Table 1).

Sulfhydryl groups may be estimated in a sample by comparison to a standard curve composed of known concentrations of a sulfhydryl-containing compound such as cysteine. Alternatively, sulfhydryl groups may be quantitated by reference to the extinction coefficient of TNB. Both methods are presented in these instructions.

Ellman's Reagent may be used for applications other than the estimation of sulfhydryls in solution. It has been used for the determination of alkylthiols by HPLC using a pre-column derivatization procedure<sup>4</sup> and to study thiols in the active site of several enzymes including thiolase,<sup>5</sup> fatty acid synthase<sup>6</sup> and mevalonate 5-diphosphate decarboxylase.<sup>7</sup>

## Procedure for Quantitating Sulfhydryl Groups Using a Cysteine Standard

### A. Material Preparation

- Reaction Buffer: 0.1M sodium phosphate, pH 8.0, containing 1mM EDTA
- Cysteine Hydrochloride Monohydrate: M.W. = 175.6, Product No. 44889 (see Related Thermo Scientific Products)
- Ellman's Reagent Solution: Dissolve 4mg Ellman's Reagent in 1mL of Reaction Buffer.

### B. Procedure

1. Prepare a set of cysteine standards by dissolving Cysteine Hydrochloride Monohydrate at the following concentrations in Reaction Buffer:

Standard	Volume of Reaction Buffer	Amount of Cysteine (M.W. = 175.6)	Final Concentration
A	100mL	26.34mg	1.5mM
B	5mL	25mL of Standard A	1.25mM
C	10mL	20mL of Standard A	1.0mM
D	15mL	15mL of Standard A	0.75mM
E	20mL	10mL of Standard A	0.5mM
F	25mL	5mL of Standard A	0.25mM
G	30mL	0ml	0.0mM (Blank)

2. Prepare a set of test tubes, each containing 50 $\mu$ L of Ellman's Reagent Solution and 2.5mL of Reaction Buffer.
3. Add 250 $\mu$ L of each standard or unknown to the separate test tubes prepared in step 2.

**Note:** For the unknown(s), make dilutions so that the 250 $\mu$ L sample applied to the assay reaction has a sulfhydryl concentration in the working range of the standard curve (0.1-1.0mM is ideal).

4. Mix and incubate at room temperature for 15 minutes.
5. Measure absorbance at 412nm.
6. Plot the values obtained for the standards to generate a standard curve. Determine the experimental sample concentrations from this curve.

**Note:** The most accurate results are obtained from the linear portion of the standard curve; i.e., the portion yielding an  $r^2$  value equal to 1.0. One or more of the high standards may exceed the linear range.

## Procedure for Quantitating Sulfhydryl Groups Based on Molar Absorptivity

### A. Material Preparation

- Reaction Buffer: 0.1M sodium phosphate, pH 8.0, containing 1mM EDTA
- Ellman's Reagent Solution: Dissolve 4mg Ellman's Reagent in 1mL of Reaction Buffer.

### B. Measure Absorbance

1. For each unknown sample to be tested, prepare a tube containing 50 $\mu$ L of Ellman's Reagent Solution and 2.5mL of Reaction Buffer.
2. Add 250 $\mu$ L of each unknown to the separate test tubes prepared in step 1. As a blank, add 250 $\mu$ L of Reaction Buffer to a separate test tube prepared in Step 1.

**Note:** For the unknown(s), make dilutions so that the 250 $\mu$ L sample applied to the assay reaction has a sulfhydryl concentration less than 1.0mM. Concentrations exceeding 1mM free sulfhydryl will result in high absorbance values and less accurate estimation of the concentration based on the extinction coefficient.

3. Mix and incubate at room temperature for 15 minutes.
4. With a spectrophotometer set to 412nm, zero the instrument on the blank and then measure absorbance of each sample.
5. Calculate the amount and concentration of sulfhydryls in the sample from the molar extinction coefficient of TNB ( $14,150\text{M}^{-1}\text{cm}^{-1}$ ), as exemplified in Section C.

### C. Example Calculation of the Free Sulfhydryl Concentration

A 250 $\mu$ L aliquot of the unknown mixed with 2.5mL of Reaction Buffer and 50 $\mu$ L of Ellman's Reagent Solution gave an absorbance of 0.879 (after subtracting the blank) using a 1cm spectrophotometric cuvette. Calculate the sulfhydryl concentration in  $\mu$ moles per mL of unknown. The reported molar absorptivity (molar extinction coefficient, which is expressed in units of  $M^{-1}cm^{-1}$ ) of TNB in this buffer system at 412nm is 14,150.<sup>2</sup> Molar absorptivity, E, is defined as follows:

$$E = \frac{A}{bc} \text{ where } A = \text{absorbance, } b = \text{path length in centimeters, } c = \text{concentration in moles/liter (=M)}$$

Solving for concentration gives the following formula:  $c = \frac{A}{bE}$

In the present example,  $A = 0.879$ ,  $b = 1cm$  and  $E = 14,150M^{-1}cm^{-1}$ . Therefore,  $c = \frac{0.879}{1(14,150)} = 6.21 \times 10^{-5} M$

This value represents the concentration of the solution in the spectrophotometric cuvette. To calculate the concentration of the unknown sample, it is necessary to account for dilution factors as follows:

The total volume of the solution being measured is

$$\begin{aligned} &2.50mL \text{ of Reaction Buffer} \\ &+ 0.25mL \text{ of Unknown Sample} \\ &+ 0.05mL \text{ of Ellman's Reagent Solution} \\ &2.80mL \text{ of solution} \end{aligned}$$

If the concentration of the assay solution is  $6.21 \times 10^{-5} M$ , then 2.80mL of that solution contains

$$2.80mL \times \frac{1L}{1000mL} \times (6.21 \times 10^{-5} \text{ moles/L}) = 1.74 \times 10^{-7} \text{ moles}$$

These  $1.74 \times 10^{-7}$  moles of sulfhydryl in the assay solution were contributed by the original 0.25mL sample. Therefore, the concentration of free sulfhydryl in the original unknown sample is

$$\frac{1.74 \times 10^{-7} \text{ moles}}{0.25mL} \times \frac{1000mL}{L} = 6.96 \times 10^{-4} M$$

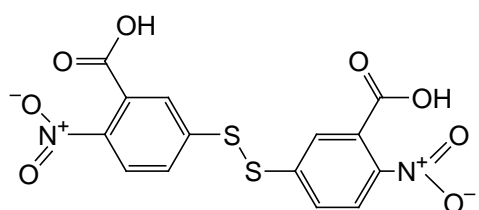
### Troubleshooting

Problem	Possible Cause	Solution
Sulfhydryl content is lower than expected	Sulfhydryls in sample are not in reduced (free) state; i.e., have become oxidized	Minimize time delay between assay of sample and its use in applications that depend on its free sulfhydryl content
		Maintain 1-5mM EDTA in sample to chelate divalent metal ions, which can oxidize sulfhydryls
		Ensure reduction of sulfhydryls using a reducing agent (see Related Thermo Scientific Products)

### Related Thermo Scientific Products

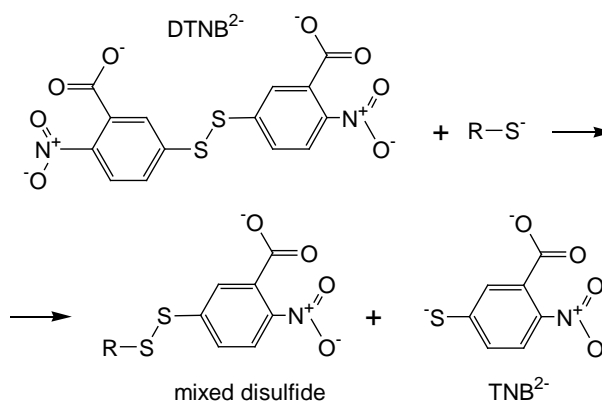
44889	Cysteine•HCl•H <sub>2</sub> O, 5g, M.W. 175.6
77712	Immobilized TCEP Disulfide Reducing Gel, 5mL
44999	SulfoLink <sup>®</sup> Immobilization Kit for Peptides, 2mL

## Additional Information



Ellman's Reagent;  
 5,5'-Dithio-bis-(2-nitrobenzoic acid)  
 $C_{14}H_8N_2O_8S_2$   
 Mol. Wt.: 396.35

**Figure 1.** Structure of Ellman's Reagent



**Figure 2.** Reduction of Ellman's Reagent

**Table 1.** Molar extinction coefficients of Ellman's Reagent in various solvents

Solvent	E at 412nm
2% SDS	12,500
0.1M phosphate, pH 8.0, 1mM EDTA	14,150
Buffered 6M guanidine hydrochloride	13,700
6M guanidine hydrochloride	13,880
8M urea	14,290

## Cited References

1. Ellman, G.L. (1959) Tissue sulfhydryl groups. *Arch Biochem Biophys* **82**:70-7.
2. Riddles, P.W., *et al.* (1983) Reassessment of Ellman's reagent. *Meth Enzymol* **91**:49-60.
3. Riddles, P.W., *et al.* (1979) Ellman's reagent: 5, 5'-dithiobis(2-nitrobenzoic acid) - a reexamination. *Anal Biochem* **94**:75-81.
4. Kuwata, K., *et al.* (1982) Liquid chromatographic determination of alkylthiols via derivatization with 5, 5'- dithiobis(2-nitrobenzoic acid). *Anal Chem* **54**:1082-7.
5. Masamune, S., *et al.* (1989) Bio-Claisen condensation catalyzed by thiolase from *Zoogloea ramigera*. Active site cysteine residues. *Chemtracts-Organic Chem* **2**:247-51.
6. Tsukamoto, Y. and Wakil, S.J. (1988) Isolation and mapping of the  $\beta$ -hydroxyacyl dehydratase activity of chicken liver fatty acid synthase. *J Biol Chem* **263**:16225-9.
7. Alvear, M., *et al.* (1989) Fractionation of chicken liver mevalonate 5-diphosphate decarboxylase by sulfhydryl-directed reagents: evidence of a functional dithiol. *Biochem Biophys Acta* **994**:7-11.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

**No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).**

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.