

Amersham Cy 5 Mono-Reactive Dye Pack

Product Specification Sheet

Introduction

Product code

PA25001

About

Reagents for the labeling of biological compounds with CyTM5 monofunctional dye.

Important

Read these instructions carefully before using the products.

Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

Storage

Store refrigerated at 2–8°C in the dark. Do not use if desiccant capsule in foil pack is either pink or green.

Expiry

See outer packaging.

Safety warnings and precautions

All chemicals should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.



NOTICE

This dye is intensely colored and very reactive. Care should be exercised when handling the dye vial to avoid staining clothing, skin and other items.

Components

- Cy5 mono-reactive dye 5 foil packs each containing sufficient dried dye to label 1 mg of protein.
- Instructions for use.

Other materials required

- **Conjugation buffer:** 0.1 M Sodium Carbonate buffer (pH 9.3).

- **Separation column:** containing a filtration gel (SephadexTM, G-25 for antibodies and G-50 for oligonucleotides).
- **Separation buffer:** phosphate-buffered saline, pH 7.2, containing 0.1% Sodium Azide.
- **Test tubes.**
- **Transfer pipettes.**
- **Glassware.**

Description

Cyanine reagents have been shown to be useful as fluorescent labels for biological compounds (1, 5). These dyes are intensely fluorescent and highly water soluble, providing significant advantages over other existing fluorophores (4).

Cy5 dye produces an intense signal in the far-red region of the spectrum. Though not recommended for visual applications, this dye is ideally suited for detection using CCD cameras, PMTs and some red-sensitive film. The Cy5 dye supplied here is a monofunctional NHS-Ester, and is provided in a dried, pre-measured form ready for the labeling of compounds containing free amino groups.

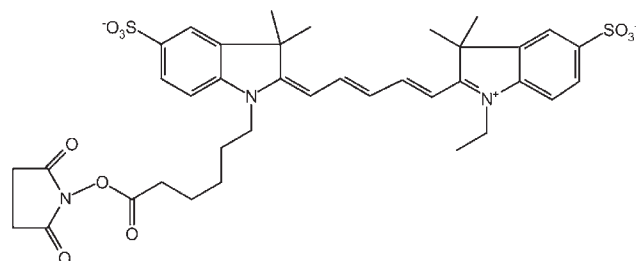


Figure 1. Cy5 monofunctional dye

Recommended procedure for use

This protocol has been designed for the preparation of Cy5-labeled IgG antibodies. It is designed to label 1 mg protein to a final molar dye/protein (D/P) ratio between 4 and 12. This assumes an average protein molecular weight of 155 000 Daltons. Other D/P ratios can be obtained by using different amounts of protein.

Note: *The following materials and procedures have been optimized for IgG antibodies. Other proteins may also be readily labeled, however, choice of buffers, separation media and technique may vary in order to produce optimal results.*

Altering the protein concentration and reaction pH will change the labeling efficiency of the reaction. Optimal labeling generally occurs at pH 9.3. Proteins have been successfully labeled with this dye at a pH as low as 7.3, however, labeling times must be significantly longer at lower pH. Higher protein concentrations usually increase labeling efficiency. Solutions of up to 10 mg/mL protein have produced good conjugation reactions.

Conjugation of dye to antibody

Antibody to be conjugated should be dissolved at 1 mg/mL in Sodium Carbonate-Sodium Bicarbonate buffer (2). Add the protein solution (1 mL) to the dye vial, cap the vial, and mix thoroughly. Care should be taken to prevent foaming of the protein solution. Incubate the reaction at room temperature for 30 minutes with additional mixing approximately every 10 minutes.

Note: Buffers containing primary amino groups such as TRIS and glycine will inhibit the conjugation reaction.

The presence of low concentrations (<2%) of biocides such as Azide or Thimerosal do not affect protein labeling.

Separation of protein from free dye

Labeled antibody can be separated from the excess, unconjugated dye by gel permeation chromatography. It is convenient to preequilibrate the column with phosphate-buffered saline and to elute the protein using the same buffer. Two blue bands should develop during elution. The faster moving band is Cy5-labeled antibody whilst the slower band is free dye. Many Cy5-labeled proteins can be stored at 2–8°C without further manipulation.

Labeled antibody can also be separated from unconjugated dye by dialysis. Dialysis does not give as efficient and rapid a separation as gel filtration. We therefore recommend that protein purification by gel filtration be used.

Estimation of final dye/protein (D/P) ratio

Dilute a portion of the labeled protein solution so that the maximum absorbance is 0.5 to 1.5 AU. Molar concentrations of dye and protein are calculated, and the ratio of these values is the average number of dye molecules coupled to each protein molecule. Molar extinction coefficients of 250 000 M⁻¹ cm⁻¹ at 650 nm for the Cy5 dye and 170 000 M⁻¹ cm⁻¹ at 280 nm for the protein are used in this example. The extinction coefficient will vary for different proteins. The calculation is corrected for the absorbance of the dye at 280 nm (approximately 5% of the absorbance at 650 nm).

$$[\text{Cy5 dye}] = (A_{650}) / 250\,000$$

$$[\text{antibody}] = (A_{280} - (0.05 \cdot (A_{650}))) / 170\,000$$

$$(D/P)_{\text{final}} = [\text{dye}] / [\text{antibody}]$$

$$(D/P)_{\text{final}} = (0.68 \cdot (A_{650})) / [A_{280} - (0.05 \cdot (A_{650}))]$$

Conjugation of dye to oligonucleotides

Modified oligonucleotides containing alkyl amino groups can be labeled with cyanine dye. Synthetic oligonucleotides must be deprotected before conjugation. Procedures that use concentrated Ammonium Hydroxide require the following pretreatment to remove all traces of Ammonia.

Concentrate the sample until it is dry (a vacuum concentrator works effectively).

Dissolve the sample in 0.25 mL of a 0.5 M Sodium Chloride solution and separate using an appropriate desalting column (Sephadex™ G-50) equilibrated with a 5.0 mM borate buffer solution adjusted to a pH of 8.0. Elute the sample with above borate buffer solution.

Concentrate the sample until it is dry. Dissolve the dry sample in a 0.1 M carbonate buffer (pH 8.5–9.0). Conjugation is carried out by adding 30 nmoles of oligonucleotide sample in approximately 0.5 mL of carbonate buffer to the dye vial. Cap the vial and mix thoroughly. Incubate the reaction at room temperature for 60 minutes with additional mixing at 15 minute intervals.

Separation of labeled oligonucleotides

Conjugated oligonucleotides can be separated from free dye using the same gel filtration procedures listed for separating conjugated antibody. A gel with a smaller exclusion size (such as Sephadex G-50) and a longer column length must be used with shorter oligonucleotides in order to ensure complete separation. Cy5-labeled oligonucleotides can be separated from unconjugated oligonucleotides using RP-HPLC. The general procedure listed in reference 3 may be optimized for the specific nucleotide sequence and HPLC configuration.

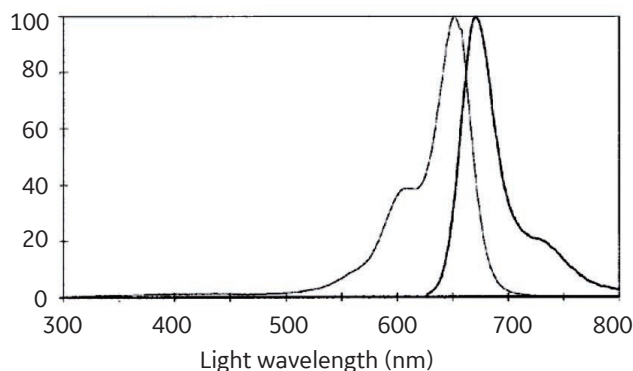


Figure 2. Cy5 dye absorption and fluorescence spectra

Cy5 monofunctional dye characteristics

Formula weight	791.99
Absorbance max	649 nm
Extinction max	250 000 M ⁻¹ cm ⁻¹
Emission max	670 nm
Quantum yield	>0.28 ¹

¹ for labeled proteins, D/P = 2

References

1. Mujumdar, R.B. *et al.*, *Bioconjugate Chemistry* **4** (2), 105-111 (1993).
2. Southwick, P.L. *et al.*, *Cytometry* **11**, 418-430 (1990).
3. Smith, L.M. *et al.*, *Nucleic Acids Research* **13**, 2399-2412 (1985).
4. Wessendorf, M.W. and Brelje, T.C., *Histochemistry* **98** (2), 81-85 (1992).
5. Yu, H. *et al.*, *Nucleic Acids Research* **22** (15), 3226-3232 (1994).

Related products

Cy3 Mono-Reactive Dye Pack	PA23001
Cy3 Maleimide Mono-Reactive Dye 5-Pack	PA23031
Cy5 Maleimide Mono-Reactive Dye 5-Pack	PA25031

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Amersham™, Cy, and Sephadex are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

All other third-party trademarks are the property of their respective owners.

The purchase of CyDye products includes a limited license to use the CyDye products for internal research and development but not for any commercial purposes. Cy and CyDye are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva. A license to use the Cy and CyDye products for commercial purposes is subject to a separate license agreement with Cytiva. Commercial use shall include:

1. Sale, lease, license or other transfer of the material or any material derived or produced from it.
2. Sale, lease, license or other grant of rights to use this material or any material derived or produced from it.
3. Use of this material to perform services for a fee for third parties, including contract research and drug screening.

If you require a commercial license to use the Cy and CyDye trademarks, please contact LSlicensing@cytiva.com.

© 2020 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

29118560 AG V:7 11/2020

