

DAPI

(4', 6-diamidino-2'-phenylindole, dihydrochloride)

62247 62248

2244.1

Number	Description
62247	DAPI, 10 mg
62248	DAPI (1 mg/ml in water), 1 ml Molecular weight: 350.25 Molar extinction coefficient: 30,600 M ⁻¹ cm ⁻¹ at 347 nm in methanol Excitation wavelength: 341 nm Emission wavelength: 452 nm CAS Number: 28718-90-3

Storage: Upon receipt store product at 2-8°C protected from light. Product is shipped at ambient temperature.

Introduction

DAPI (4', 6-diamidino-2'-phenylindole, dihydrochloride) is a dye that fluoresces blue (455 nm) when bound to double-stranded DNA and excited by exposure to 345 nm light. DAPI binds selectively to the minor groove of adenine-thymine (A-T) regions of DNA, where its fluorescence is approximately 20-fold greater than in the nonbound state.¹⁻³ These properties make DAPI useful for assaying DNA in solution,⁴ diagnosing mycoplasma infection of cell cultures,⁵ measuring nuclear content and sorting chromosomes in flow cytometry,⁶ assessing apoptosis,⁷ and detecting nuclei and organellar DNA in immunofluorescent and *in situ* hybridization procedures.^{2,8} DAPI is also used as a replacement for ethidium bromide to stain double-stranded DNA in agarose gels.^{5,9} Because DAPI is cell permeable and fluoresces blue, it is commonly used to counterstain nuclei in histochemical methods when red-fluorescent antibodies have been used to detect specific targets.⁸ Reports also indicate that DAPI binds to polyphosphates and other polyanions,¹⁰ dextran sulfate¹¹ and SDS.¹²

General Procedure for Immunofluorescent Staining

A. Reagent Preparation

PBS (Wash Buffer)	Modified Dulbecco's PBS (Product No. 28374): 8 mM sodium phosphate, 2 mM potassium phosphate, 140 mM sodium chloride, 10 mM potassium chloride; pH 7.4.
DAPI Stock Solution	Dissolve DAPI in ultrapure water to 1 mg/ml. Stock solution is stable for several months and repeated use if stored protected from light at -20°C.
DAPI Working Solution	Dilute the DAPI Stock Solution 1:1,000 in ultrapure water or PBS (1 µg/ml DAPI). Filter the Working Solution to remove dye aggregates that can result in punctate signal.

B. Procedure

1. Follow standard procedures to fix sample and then probe with specific fluorescent-labeled antibodies.
2. Thoroughly wash sample with PBS to remove nonbound probe.
3. Add a sufficient volume of DAPI Working Solution to completely cover the sample. Place aluminum foil over the sample to protect it from light and incubate at room temperature for 2-10 minutes. If resulting final signal bleeds through to other fluorescent channels, decrease incubation time in next experiment.
4. Wash sample thoroughly with PBS to remove excess DAPI.
5. Mount sample with an appropriate medium and detect according to standard protocols.

Procedure for Assaying DNA

DAPI can be used to quantitate DNA in solution. The method is relatively insensitive to pH 5-10 but is sensitive to changes in temperature and ionic strength, as well as to fluorescence quenching by divalent or heavy metal cations.⁴ The fluorescence is not linear over broad concentration ranges of DNA; therefore, use an internal standard each time the assay is performed.

A. Reagent Preparation

Assay Buffer	0.1 M NaCl, 10 mM EDTA, 10 mM Tris; pH 7.0.
DAPI Stock Solution	Dissolve DAPI in ultrapure water to 1 mg/ml. Stock solution is stable for several months and repeated use if stored protected from light at -20°C.
DAPI Working Solution	0.1 µg/ml DAPI in Assay Buffer.
DNA Standards	Dilute known amount of calf thymus DNA with Assay Buffer to make a series of DNA standards with concentrations ranging from 0 to 5 µg/ml (0 to 250 ng/50 µl). Prepare replicates of each dilution so error statistics can be calculated.

B. Assay

1. Add 50 µl of each unknown sample or DNA Standard to a disposable fluorometer cuvette.
2. Add 1.5 ml DAPI Working Solution to each cuvette.
3. Cover cuvettes with foil and incubate at room temperature for 10 minutes.
4. Measure fluorescence of each solution.
5. Prepare a standard curve by plotting the mean values of the standards plotted against their concentrations.
6. Determine concentration of each unknown sample based on its fluorescence measurement relative to the standard curve.

Related Thermo Scientific Products

Visit our web site for information about other fluorescent protein labeling reagents, including fluorescein (Product No. 46409, 46410 and 53029), rhodamine (No. 46406 and 53031), and amine-reactive or sulfhydryl-reactive Thermo Scientific DyLight Fluors. We also offer a variety of labeled secondary antibodies.

References

1. Morikawa, K. and Yanagida, J. (1981). Visualization of individual DNA molecules in solution by light microscopy: DAPI staining method. *J. Biochem. (Tokyo)* **89**:693-6.
2. Lawrence, M.E. and Possingham, J.V. (1986). Direct measurement of femtogram amounts of DNA in cells and chloroplasts by quantitative microspectrofluorometry. *J. Histochem. Cytochem.* **34**:761-8.
3. Kubista, M., *et al.* (1987). Characterization of interaction between DNA and 4', 6-diamidino-2-phenylindole by optical spectroscopy. *Biochemistry* **26**:4545-53.
4. Brunk, C.F., *et al.* (1979). Assay for nanogram quantities of DNA in cellular homogenates. *Anal. Biochem.* **92**:497-500.
5. Russell, W.C., *et al.* (1975). A simple cytochemical technique for demonstration of DNA in cells infected with mycoplasmas and viruses. *Nature* **253**:461-2.
6. Hammarton, T.C., *et al.* (2003). Stage-specific differences in cell cycle control in *Trypanosoma brucei* revealed by RNA interference of a mitotic cyclin. *J. Biol. Chem.* **278**(25):22877-86.
7. Lai, J., *et al.* (2003). Loss of HSulf-1 up-regulates heparin-binding growth factor signaling in cancer. *J. Biol. Chem.* **278**(25):23107-17.
8. Soto, P., *et al.* (2003). SMAD2 and SMAD7 Involvement in the post-translational regulation of Muc4 via the transforming growth factor-β and interferon-γ pathways in rat mammary epithelial cells. *J. Biol. Chem.* **278**(22):20338-44.
9. Nairn, R.S., *et al.* (1982). Comparison of ethidium bromide and 4', 6'-diamidino-2-phenylindole as quantitative fluorescent stains for DNA in agarose gels. *J. Biochem. Biophys. Meth.* **6**:95-103.
10. Tijssen, J.P.F., *et al.* (1982). Localization of polyphosphates in *Saccharomyces fragilis*, as revealed by 4', 6-diamidino-2-phenylindole fluorescence. *Biochem. Biophys. Acta* **721**:394-8.
11. Allan, R.A. and Miller, J.J. (1980). Influence of S-adenosylmethionine on 4', 6-diamidino-2-phenylindole dihydrochloride (DAPI)-induced fluorescence of polyphosphate in the yeast vacuole. *Can. J. Micro.* **26**:912-20.
12. Kapuscinski, J. and Skoczylas, B. (1978). Fluorescent complexes of DNA with DAPI (4', 6-diamidine-2-phenyl indole dihydrochloride) or DCI (4', 6-dicarboxamide-2-phenyl indole). *Nucl. Acids Res.* **5**:3775-99.

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 versions of product instructions are available at www.thermo.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2010 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.