

Singlet Oxygen Sensor Green Reagent

S36002 Singlet Oxygen Sensor Green Reagent *special packaging*

Quick Facts

Storage upon receipt:

- $\leq -20^{\circ}\text{C}$
- Protect from light
- Desiccate

Ex/Em of reaction product: 504/525 nm

Introduction

The Singlet Oxygen Sensor Green reagent is highly selective for $^1\text{O}_2$; unlike other available fluorescent and chemiluminescent singlet oxygen detection reagents, it does not show any appreciable response to hydroxyl radical ($\bullet\text{OH}$) or superoxide ($\bullet\text{O}_2^-$) (Figure 1). This new singlet oxygen indicator initially exhibits weak blue fluorescence, with excitation peaks at 372 and 393 nm and emission peaks at 395 and 416 nm. In the presence of singlet oxygen, it emits a green fluorescence similar to that of fluorescein (excitation/emission maxima $\sim 504/525$ nm). The Singlet Oxygen Sensor Green reagent is supplied as a cell-impermeant derivative.

Materials

Contents

Singlet Oxygen Sensor Green reagent is provided specially packaged in sets of 10 vials, each containing 100 μg .

Storage and Handling

Upon receipt, store the Singlet Oxygen Sensor Green reagent desiccated and protected from light at $\leq -20^{\circ}\text{C}$ until required for use. Prepare stock solutions in methanol; dissolve the contents of one 100 μg vial in 33 μL of methanol to make a stock solution of ~ 5 mM. Prepare working solutions of this reagent immediately before use, and discard any excess diluted reagent at the end of the work session.

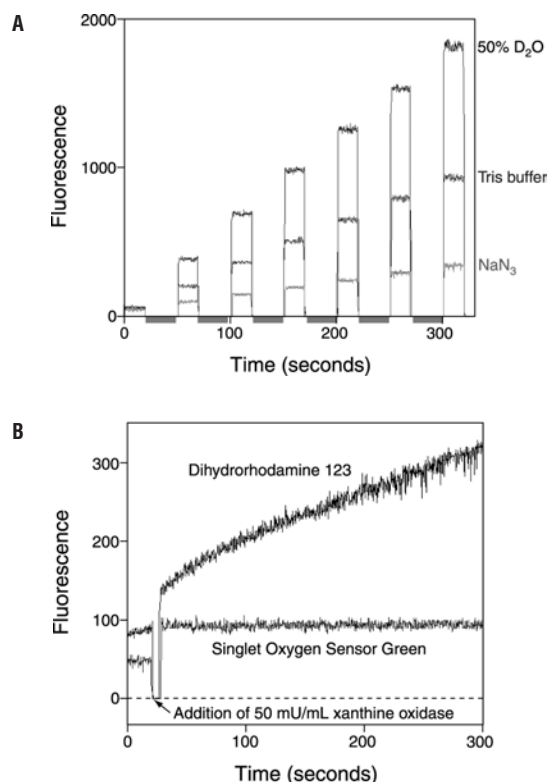


Figure 1. Fluorescence response and specificity of Singlet Oxygen Sensor Green reagent to $^1\text{O}_2$. (A) Fluorescence measurements were made in a spectrofluorometer using excitation/emission of 488/525 nm for solutions containing: 1 μM Singlet Oxygen Sensor Green reagent and 10 μM methylene blue in 100 mM pH 7.5 Tris buffer alone; the singlet oxygen scavenger sodium azide (NaN_3); or 50% D_2O , which increases the lifetime of $^1\text{O}_2$. Measurements were made for 20-second periods, with 30-second intervals (indicated by grey bars) between each measurement. During the 30-second intervals, the samples were exposed to laser radiation (630–680 nm, <5 mW), resulting in methylene blue-photosensitized generation of $^1\text{O}_2$. (B) Fluorescence measurements were made in a spectrofluorometer using excitation/emission of 488/525 nm for solutions of 50 mM pH 7 Tris buffer with 1 mM xanthine containing either 1 μM Singlet Oxygen Sensor Green reagent or dihydrorhodamine 123. After ~ 20 seconds, 50 mU/mL of xanthine oxidase (XO) was added. XO catalyzes the oxidation of xanthine, producing uric acid and superoxide. Superoxide can spontaneously degrade to H_2O_2 .

Application

The Singlet Oxygen Sensor Green reagent should be useful for detecting $^1\text{O}_2$ in aqueous solutions. It can also potentially be

employed to assess the efficacy of free radical scavengers, which are frequently used to improve the flavor and nutritional quality of foods. Note that the Singlet Oxygen Sensor Green reagent can also become activated at alkaline pH or in the presence of certain solvents, including (but not limited to) acetonitrile, DMSO, DMF, and acetone. Its fluorescent product may also degrade over time in some solutions. However, with proper controls, the intensity of the green-fluorescent signal can be correlated with $^1\text{O}_2$

concentration, without significant interference from other reactive oxygen species.

Solution Assays

Singlet Oxygen Sensor Green reagent is intended for use in aqueous environments. The optimal dilution buffer and working concentration should be determined empirically; a suggested starting concentration range is 1–10 μM .

Reference

1. Free Radic Biol Med 27, 873-881 (1999).

Product List *Current prices may be obtained from our Web site or from our Customer Service Department.*

Cat #	Product Name	Unit Size
S36002	Singlet Oxygen Sensor Green *special packaging*	10 x 100 μg

Contact Information

Further information on Molecular Probes' products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Leiden, the Netherlands. All others should contact our Technical Assistance Department in Eugene, Oregon.

Please visit our Web site — www.probes.com — for the most up-to-date information.

Molecular Probes, Inc.

29851 Willow Creek Road, Eugene, OR 97402
Phone: (541) 465-8300 • Fax: (541) 335-0504

Customer Service: 6:00 am to 4:30 pm (Pacific Time)
Phone: (541) 335-0338 • Fax: (541) 335-0305 • order@probes.com

Toll-Free Ordering for USA and Canada:
Order Phone: (800) 438-2209 • Order Fax: (800) 438-0228

Technical Assistance: 8:00 am to 4:00 pm (Pacific Time)
Phone: (541) 335-0353 • Toll-Free (800) 438-2209
Fax: (541) 335-0238 • tech@probes.com

Molecular Probes Europe BV

Poortgebouw, Rijsburgerweg 10
2333 AA Leiden, The Netherlands
Phone: +31-71-5233378 • Fax: +31-71-5233419

Customer Service: 9:00 to 16:30 (Central European Time)
Phone: +31-71-5236850 • Fax: +31-71-5233419
eurorder@probes.nl

Technical Assistance: 9:00 to 16:30 (Central European Time)
Phone: +31-71-5233431 • Fax: +31-71-5241883
eurotech@probes.nl

Molecular Probes' products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Several of Molecular Probes' products and product applications are covered by U.S. and foreign patents and patents pending. Our products are not available for resale or other commercial uses without a specific agreement from Molecular Probes, Inc. We welcome inquiries about licensing the use of our dyes, trademarks or technologies. Please submit inquiries by e-mail to busdev@probes.com. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2004, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.