Probenecid, water soluble

 Table 1. Contents and storage information.

Material	Amount	Storage	Stability
Probenecid, water soluble	10 × 77 mg	 ≤25°C Desiccate 	When stored as directed the product is stable for 6 months

Introduction

Probenecid is commonly used to inhibit organic-anion transporters located in the cell membrane. Such transporters can extrude dyes and indicators and thus contribute to poor loading or a high background signal in assays based on retention of the dyes or indicators inside cells. For example, fluo-4, AM is a Ca^{2+} indicator widely used for in-cell measurement of agonist-stimulated and antagonist-inhibited calcium signaling. The acetoxymethyl ester (AM) dye precursor is uncharged, cell permeant, and nonfluorescent. Inside the cell, nonspecific esterases cleave the AM blocking groups, generating the charged, active form of fluo-4 which fluoresces upon Ca^{2+} binding. Although the charged dye molecules leak out of cells far more slowly than the uncharged precursor molecules, the anion transporters promote the leakage of the fluorescent dye over time. The use of probenecid to block the efflux of intracellular dyes was first reported by Di Virgilio et al.¹

Wash steps or quencher dyes may be incorporated into fluorescent assays to minimize baseline fluorescence. However, washing introduces an extra step that is undesirable for high-throughput applications, and may also risk loss of non-adherent cells. Quencher dyes, while offering the advantage of homogeneous (one-step, mix-and-read) assays, may interact negatively with some receptor systems of interest.² The use of probenecid to suppress efflux of fluorescent dyes is a favorable method for reducing baseline fluorescence. Our water-soluble probenecid, available as a stand-alone product or supplied with our Fluo-4 NW (no-wash) Calcium Assay Kits, has the advantages of being easy to dissolve in buffer and safer to use than the commonly used free acid form, which requires caustic 1 M NaOH to dissolve it. It is also slightly more effective than the free acid form at equimolar concentrations, probably due to its better solubility (Figure 1).

Guidelines for Use

Using Probenecid

Store the dry, solid probenecid desiccated at $\leq 25^{\circ}$ C. Dissolve the contents of one vial in 1 mL of buffer, such as HBSS, and use the solution the same day or store it at $\leq -20^{\circ}$ C for up to 6 months. This stock solution of probenecid is at a concentration of 250 mM. Typical working concentrations in cell-based assays are 1–2.5 mM.

References

1. Cell Calcium 11, 57 (1990). 2. Biotechniques 34, 164 (2003).



Figure 1. Inhibition of dye extrusion using 2.5 mM probenecid. CHO M1 cells were incubated in dye loading solution for 60 minutes at 37°C, which accentuates the probenecid effect. Calcium release was triggered in CHO M1 cells (M1WT3 cell line, ATCC #CRL-1985) using 20 nM carbachol. The averages of four measurements (without baseline subtraction) in a Fluo-4 NW assay were plotted.

Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name U	nit Size
P36400	Probenecid, water soluble	< 77 mg
P10020	PowerLoad™ Concentrate, 100X	5 mL
F36205	Fluo-4 NW Calcium Assay Kit (high-throughput) *for 100 microplates*	1 kit
F36206	Fluo-4 NW Calcium Assay Kit (starter pack with buffer) *for 10 microplates*	1 kit

Contact 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 probesorder@invitrogen.com

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

Technical Service:

8:00 am to 4:00 pm (Pacific Time) Phone: (541) 335-0353 Toll-Free (800) 438-2209 Fax: (541) 335-0238 probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd. 3 Fountain Drive Inchinnan Business Park Paisley PA4 9RF, UK Phone: +44 (0) 141 814 6100 Fax: +44 (0) 141 814 6260 Email: euroinfo@invitrogen.com Technical Services: eurotech@invitrogen.com 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 Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

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.

Limited Use Label License No. 223: Labeling and Detection Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation [®] are registered with the U.S. Patent and Trademark Office.

Copyright 2008, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.