INSTRUCTIONS p-Bromophenacyl-8TM Reagent



1521.0

TS-48891

Description

Number TS-48891

p-Bromophenacyl-8[™] Reagent, 10 ml Hypo-Vial[™] Storage Vial, contains 0.1 mmol/ml *p*bromophenacyl bromide and 0.005 mmol/ml crown ether in acetonitrile Molecular Weight: 277.94

λmax: 260 nm

Storage: Upon receipt store product at room temperature.

Introduction

p-Bromophenacyl-8TM Reagent is an excellent derivatization reagent that reacts with carboxylic acids. This reagent is used for preparing phenacyl esters, which are used to separate many saturated and unsaturated fatty acids¹⁻² for UV detection in HPLC applications with low nanomole sensitivity. *p*-Bromophenacyl-8TM Reagent is also useful for studying acid composition of bacterial cell walls.³⁻⁴

Procedure for Preparing Phenacyl Esters

1. Dissolve ~10 mg acid in MeOH in a 5.0 ml Reacti-Vial[™] Small Reaction Vial (Product No. TS-13223) fitted with a magnetic stirrer (Product No. TS-16000). Neutralize to the phenolphthalein endpoint with 85% KOH in MeOH.

Note: If the formation of potassium salts is undesirable, instead of using KOH, neutralize by adding KHCO₃ at 3-5 times the molar equivalent of the free acid.

- 2. Evaporate the MeOH with N₂.
- 3. Add 1.0 ml of *p*-Bromophenacyl-8TM Reagent and 2.0 ml dry acetonitrile to the reaction vial.
- 4. Heat reaction at 80°C with stirring for 30 minutes.
- 5. Remove reaction vial from the heating block and allow it to cool.
- 6. Analyze solution by HPLC. Use a reverse phase C18 column with acetonitrile and water as the mobile phase.

Cited References

- 1. Ahmed, M. S., *et al.* (1980). Use of *p*-bromophenacyl bromide to enhance ultraviolet detection of water-soluble organic acids (steviolbioside and rebaudioside B) in high-performance liquid chromatographic analysis. *J. Chromatogr.* **192:**387-93.
- 2. Borch, R. F. (1975). Separation of long chain fatty acids as phenacyl esters by high pressure liquid chromatography. Anal. Chem. 47(14):2437-9.
- 3. Manzoor, S. E., *et al.* (1999). Reduced glutaraldehyde susceptibility in *Mycobacterium chelonae* associated with altered cell wall polysaccharides. J. Antimicrob. Chemother. **43**:759-765.
- 4. Sokolovská, I., *et al.* (2003). Carbon source-induced modifications in the mycolic acid content and cell wall permeability of *Rhodococcus erythropolis* E1. *Appl. Environ. Microbiol.* **69**(**12**):7019-27.

General Reference

Knapp, D. R. (1979). Handbook of analytical derivatization reactions. John Wiley and Sons, New York; Chapter 3.

Current versions of product instructions are available at <u>www.thermo.com/columns</u>. © 2008 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.

Warranty: Pierce Biotechnology products are warranted to meet stated product specifications and to conform to label descriptions when stored and used properly. Unless otherwise stated, this warranty is limited to one year from date of sale when used according to product instructions. Pierce Biotechnology's sole liability for the product is limited to replacement of the product or refund of the purchase price. Unless otherwise expressly authorized in writing by Pierce Biotechnology, Pierce products are supplied for Research Use Only and are intended to be used by a technically qualified individual. Pierce Biotechnology's quality system is certified to ISO 9001. Pierce Biotechnology products are not produced in accordance with FDA's current Good Manufacturing Practices. Pierce Biotechnology strives for 100% customer satisfaction. If you are not satisfied with the performance of a Pierce Biotechnology product, please contact Pierce Biotechnology or your local distributor.