

# *p*-Bromophenacyl-8<sup>TM</sup> Reagent

48891

1521.0

**Number****Description**

48891

***p*-Bromophenacyl-8<sup>TM</sup> Reagent**, 10 ml Hypo-Vial<sup>TM</sup> Storage Vial, contains 0.1 mmol/ml *p*-bromophenacyl bromide and 0.005 mmol/ml crown ether in acetonitrile

Molecular Weight: 277.94

$\lambda_{\text{max}}$ : 260 nm

**Storage:** Upon receipt store product at room temperature.

**Introduction**

*p*-Bromophenacyl-8<sup>TM</sup> 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 acids<sup>1-2</sup> for UV detection in HPLC applications with low nanomole sensitivity. *p*-Bromophenacyl-8<sup>TM</sup> Reagent is also useful for studying acid composition of bacterial cell walls.<sup>3-4</sup>

**Procedure for Preparing Phenacyl Esters**

1. Dissolve ~10 mg acid in MeOH in a 5.0 ml Reacti-Vial<sup>TM</sup> Small Reaction Vial (Product No. 13223) fitted with a magnetic stirrer (Product No. 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 KHCO<sub>3</sub> at 3-5 times the molar equivalent of the free acid.

2. Evaporate the MeOH with N<sub>2</sub>.
3. Add 1.0 ml of *p*-Bromophenacyl-8<sup>TM</sup> 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 C<sub>18</sub> 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.

---

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](http://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.