Enhancement of Detection Sensitivity of Tobramycin Using Pre-column Derivatization

Varian Application Note Number 5

Fran Lai Varian Chromatography Systems

Key Words: Pharmaceuticals, Tobramycin, OPA, AutoMix, Detection Enhancement, Fluorescence, Pre-column Derivatization, Varian 9100 AutoSampler

Introduction

Reversed Phase HPLC and UV detection are very popular methods for the analysis of pharmaceutical compounds. This works well for nonpolar and UV-absorbent compounds, but for compounds that are rather polar and lack UVabsorptivity, the analysis becomes a major challenge.

Tobramycin, an aminoglycoside antibiotic, is a good example of a polar pharmaceutical compound with low UV-absorptivity. To overcome these problems, a method has been developed using pre-column derivatization with orthophthalaldehyde (OPA).

Procedure

The derivatization process is easily automated using the sample preparation features on the 9100 AutoSampler.

- 1. Prepare Tobramycin sample in acetonitrile:water/60:40.
- 2. Prepare OPA reagent:

OPA 5 mg/mL

2-mercaptoethanol (2-ME) 1% v/v in borate buffer pH 10.4, 0.4M

Derivatization Using Varian 9100 AutoSampler

- 1. Transfer and mix OPA reagent with Tobramycin sample (1:1 v/v) (Figure 1).
- 2. Wait for programmed reaction time.
- 3. Inject onto the HPLC.

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to www.agilent.com/chem.

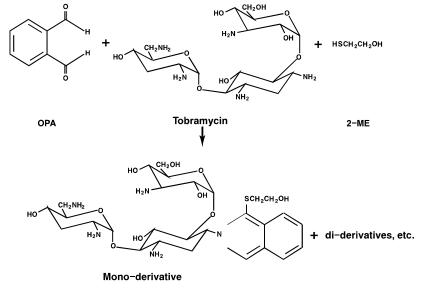
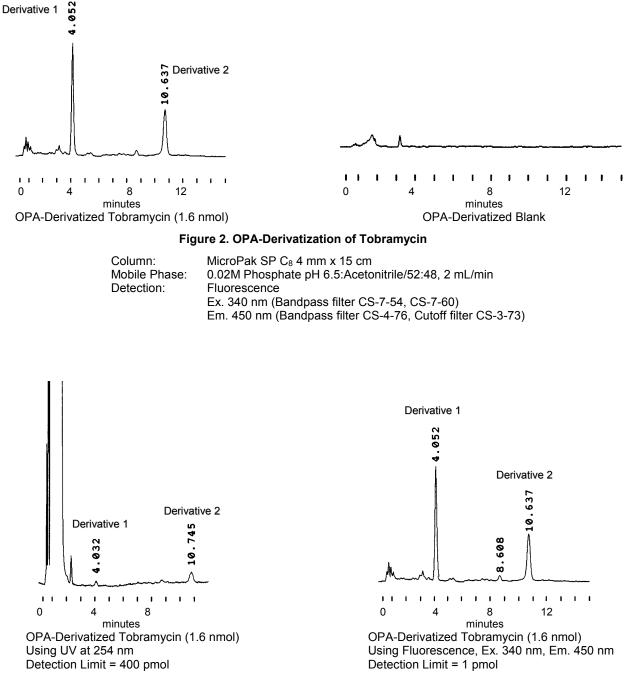


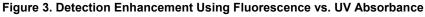
Figure 1. Analysis of Tobramycin Using Pre-column Derivatization (OPA)

Results

- 1. Two derivatives are obtained. Due to multiple primary amino sites on the compound, multiple derivatives may be expected. The ratio of Derivative 2/Derivative 1 increases with reaction time and with increase in acetonitrile in the sample solvent.
- 2. The optimum reaction time is 30 minutes, and the optimum solvent composition is 60% acetonitrile. There is no interference from the blank. (Figure 2)
- 3. Using fluorescence, the detection limit is about 400 times better than when using UV absorbance. (Figure 3)
- 4. Linearity was tested up to 1.2 nmoles and found to be linear with a correlation coefficient of 0.998.







References

- Sten-Erik Bäck, Chemical Assay, Involving Liquid Chromatography, for Aminoglycoside Antibiotics in Serum, Clinical Chemistry, Vol 25, No. 7, p. 1222-1225, 1979.
- 2. Robert Cunico, Automatic Precolumn Derivatization of Amino Acids with Orthophthalaldehyde, Varian LC at Work #161, 1985.

