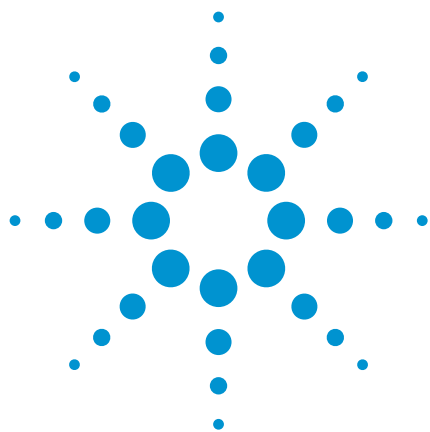




Monam in sewage water

Analysis of pesticides

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography using an Agilent CP-Sil 5 CB column isolates methyl isothiocyanate, an indicator of the soil fumigant Monam, in water from a sewage treatment plant in eight minutes.



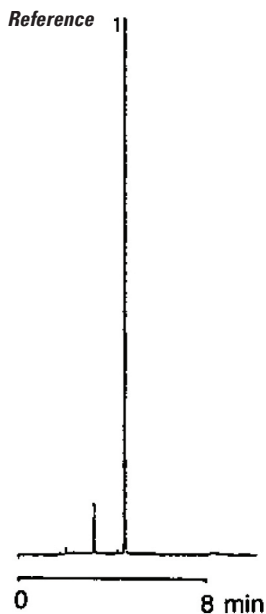
Agilent Technologies

Conditions

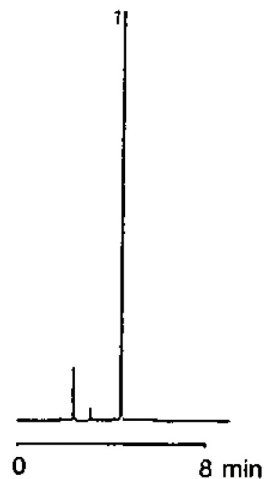
Technique	: GC-PTI
Column	: Agilent CP-Sil 5 CB, 0.32 mm x 25 m fused silica WCOT CP-Sil 5 CB (1.2 µm) (Part no. CP7760)
Temperature	: 80 °C
Carrier Gas	: N ₂ , 50 kPa (0.5 bar, 7 psi)
Injector	: PTI cryofocussing : -80 °C precool time : 2 min purge time : 10 min purge flow : 10 mL/min condenser temp. : -15 °C injection temp. : 200 °C injection time : 5 min
Detector	: FID, 8 x 10 ⁻¹² Afs
Concentration Level	: reference: 0.72 µg/L (0.72 ppb) sewage water: 0.5 µg/L (0.50 ppb)
Sample Size	: reference: 5 mL (diluted 1:2 x 10 ⁶) sewage water: 5 mL (diluted 1:10 ³)
Courtesy	: R. Dilling, H. de Bie and G. Sibon, Zuiveringsschap Drenthe, Assen, The Netherlands

Peak identification

1. methyl isothiocyanate



Sewage water sample



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A00298



Agilent Technologies