Automatic Precolumn Derivatization for the HPLC Determination of Aliphatic Amines in Air

Huang Xiongfeng,¹ Xu Qun,¹ and Jeffrey Rohrer²
¹Thermo Fisher Scientific, Shanghai, People's Republic of China; ²Thermo Fisher Scientific, Sunnyvale, CA, USA

Key Words

Volatile Aliphatic Amines, Environmental Analysis, Air Pollutant, Acclaim PolarAdvantage II Column, Fluorescence Detection

Introduction

Aliphatic amines such as methylamine, ethylamine, diethylamine, propylamine, and butylamine are considered important air pollutants due to their odorous and toxic properties. They are found in the air as a result of their industrial commercial applications and their widespread uses as intermediates in the chemical and pharmaceutical industries. In addition to their toxicity, these amines have been recognized as potential precursors of carcinogenic nitrosamines in the presence of nitrous acid or other nitrosating agents in polluted air. Therefore, determination of amines at low concentrations in the atmosphere is justified by growing concern about the public health hazards that these compounds present.

The techniques most widely used for the determination of aliphatic amines are gas chromatography (GC), ^{1,3-6} high-performance liquid chromatography (HPLC),^{2,7-13} ion chromatography,¹⁴⁻²¹ and spectrophotometry.^{22,23} Because these amines are subject to adsorption and decomposition, it is preferable to derivatize them to a stable product prior to analysis.

Precolumn chemical derivatization in solution has been accepted as an effective modification technique prior to HPLC analysis. This technique can provide highly sensitive and selective detection of these amines by bonding a chromophore or fluorophore that results in products with strong UV absorption and/or fluorescence emission. This approach can also improve retention by reducing the polarity of these compounds.

Many derivatization reagents for primary and secondary aliphatic amines prior to HPLC analysis have been reported, such as o-phthalaldehyde (OPA), 2, 4-dinitrofluorobenzene (DNFB), 9-fluorenylmethyl chloroformate (FMOC-Cl), dinitrochlorobenzene (DNCB), 5-dimethylaminonaphthalene-1-sulfonyl chloride (DNS-Cl), and phenylisothiocyanate (PITC).24 FMOC-Cl and OPA are often used due to the rapid formation of stable derivatives under mild conditions—preferably at room temperature and in aqueous phase—and have been applied to the determination of amino acids with automatic precolumn derivatization.²⁵ Automatic precolumn derivatization eliminates tedious manual procedures, reduces error, and thereby increases method reproducibility. This technique is performed using the autosampler and controlled by the chromatography data system software.

Goal

To develop an efficient HPLC method using automatic precolumn derivatization and fluorescence detection for the determination of ethanolamine, methylamine, ethylamine, diethylamine, *n*-propylamine, *n*-butylamine, and *n*-hexylamine in air samples



Equipment

- Thermo Scientific Dionex UltiMate 3000 HPLC system, including:
 - DGP-3600RS Dual Ternary Rapid Separation Pump System with SRD-3600 Integrated Solvent and Degasser Rack
 - WPS-3000TRS Rapid Separation Wellplate Sampler, Thermostatted, with 250 µL sample loop
 - TCC-3000RS or TCC-3000SD Thermostatted Column Compartment
 - FLD-3400RS Rapid Separation Fluorescence Detector
- Thermo Scientific Dionex Chromeleon Chromatography Data System (CDS) software version 6.80, SR9 or higher
- Thermo Scientific Orion 2-Star Benchtop pH Meter
- SK3200 KUDOS LHC Series Dual Frequency Ultrasonic Cleaner, KUDOS, Shanghai, China

Reagents and Standards

- Deionized (DI) water, 18.2 M -cm resistivity
- Methanol (CH₃OH), HPLC Grade (Fisher Chemical P/N AC610090040)
- Acetonitrile (CH₃CN), HPLC Grade (Fisher Chemical P/N AC610010040)
- Phosphoric Acid (H₃PO₄), Analytical Grade, SCRC, China
- Sodium Dihydrogen Phosphate (NaH₂PO₄), Analytical Grade, SCRC, China
- Sodium Hydroxide (NaOH), Analytical Grade, SCRC, China
- 9-Fluorenylmethyl Chloroformate (FMOC-Cl), ≥99.0% (HPLC), Sigma-Aldrich®
- *o*-Phthalaldehyde (OPA) ≥99.0% (HPLC), Sigma-Aldrich
- 3-Mercaptopropionic Acid ≥99.0%, Alfa, Tianjin, China
- Sodium Borate (Na₂B₄O₇·10H₂O), Analytical Grade, SCRC, China
- Ethanolamine, Methylamine, Ethylamine, Diethylamine, *n*-Propylamine, *n*-Butylamine, and *n*-Hexylamine, Analytical Grade, SCRC, China

Solutions for Derivatization

Borate buffer (400 mM, pH 10.2): dissolve 15.26 g of $Na_2B_4O_7\cdot 10H_2O$ in 50 mL DI water, then dilute to 100 mL with DI water. Adjust to pH 10.2 with 1 M NaOH.

Borate buffer (20 mM, pH 10.2): dilute 5 mL borate buffer (400 mM) to 100 mL with DI water, then adjust to pH 10.2 with 1 M NaOH.

Phosphate buffer (50 mM, pH 3.0, used as injection diluent): dissolve 1.53 g of NaH_2PO_4 in 100 mL DI water, then dilute to 250 mL with DI water. Adjust to pH 3.0 with 50% H_3PO_4 .

OPA stock solution (50 mg/mL): dissolve 50 mg of OPA in 1 mL methanol.

OPA solution (75 mg/L): dilute 75 μ L OPA stock solution (50 mg/mL) to 50 mL with methanol.

OPA reagent: combine 0.8 mL 20 mM borate buffer + 0.2 mL 75 mg/L OPA + 2 µL 3-mercaptopropionic acid.

FMOC reagent (2.5 mg/mL): dissolve 50 mg of FMOC in 20 mL acetonitrile.

Note: The OPA and FMOC solutions and reagents are sensitive to oxidation and light; therefore, for optimum sensitivity they should be freshly prepared in reagent reservoirs that have been rinsed with methanol.

Preparation of Mobile Phase

Phosphate buffer (50 mM, pH 8.5): dissolve 3.0 g of NaH_2PO_4 in 500 mL water, then adjust to pH 8.5 with 1 M NaOH.

Working Standard Solutions for Calibration

Prepare each amine standard by accurately weighing 25 mg of a standard and diluting to 250 mL in a volumetric flask with water/methanol (1:1, v/v) solution. The concentration of the standard is 100 mg/L (stock standard solution). Prepare four working standard solutions for the calibration with 0.1, 0.5, 1.0, 2.0, and 5.0 μ g/mL concentrations by adding the proper amount of stock standard solution and diluting with a water/ methanol (1:1, v/v) solution.

Sample Preparation

Air Sample Extraction

Procedures for the extraction are based on those specified in a publication.²⁶ In this study, particulate matter (≤2.5 μm) in the atmosphere was collected on a quartz membrane (Ø 90 mm, Waterman, UK) for 48 h at a flow rate of 77.49 L/min using a TSP/PM2.5 sampler (Dike Mechanical & Electrical Technology Ltd. Co., Beijing, China) that was located at the top of Building No. 4, Fudan University, Shanghai, China. The compounds of interest were then extracted with 10 mL of water/methanol (1:1, v/v) solution in an ultrasonic bath for 40 min and cooled to room temperature. The extract was filtered through a 0.45 μm membrane (Millex™-LHR) filter prior to derivatization.

Sample Derivatization

The derivatization reactions for primary and secondary aliphatic amines using OPA and FMOC are shown in Figure 1.24 Place the derivatization reagents, borate buffer (400 mM), OPA reagent (0.8 mL 20 mM borate buffer + 0.2 mL 75 mg/L OPA + 2 μL 3-mercaptopropionic acid), FMOC reagent (2.5 mg/mL), and injection dilution buffer (50 mM phosphate buffer, pH 3.0) each in different vials in the autosampler. Automatically execute the derivatization by drawing the derivatization reagents and sample solutions into the sample loop. First, draw reagents and sample solutions into the sample loop in the order of borate buffer (50 µL), OPA reagent (2.5 µL), and sample solution (1 µL); then move the syringe several times to mix the solution. This part of the procedure is for the derivatization of the primary amines. Second, draw FMOC reagent (1 µL) into the sample loop and mix by moving the syringe several times to derivatize the secondary amines.

After derivatization, the derivatized sample solutions are automatically flushed from the sample loop into the analytical flow for amine separations and detection. The whole process is controlled by Chromeleon™ CDS software using the program as shown in Table 1.

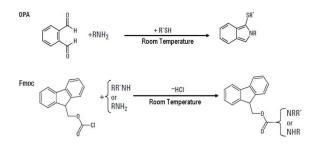


Figure 1. Derivatization reactions of primary and secondary aliphatic amines using OPA and FMOC.

Conditions				
Column:	Thermo Scientific Acclaim PolarAdvantage II (PA2), 3 μ m Analytical (3.0 \times 150 mm)			
Mobile Phase:	A: 50 mM phosphate buffer, pH 8.5 B: CH ₃ CN/CH ₃ OH (1:1, v/v)			
Gradient:	B: 0–4 min, 35%; 4–7 min, 35-55%; 7–15 min, 55%			
Flow Rate:	0.6 mL/min			
Temperature:	40 °C			
Detection:	Fluorescence, Ex: 230 nm, Em. 450 nm			

Step	Program	Note		
	DrawSpeed = 10.000 [μL/s]			
	DrawDelay = 3000 [ms]			
	DispSpeed =10.000 [μ L/s]			
	DispenseDelay = 0 [ms]			
	WasteSpeed =10.000 [µL/s]			
	SampleHeight = 2.000 [mm]			
	InjectWash = Both	D 100 100 100 100 100 100 100 100 100 10		
	WashVolume = $250.000 [\mu L]$	ReagentAVial: 400 mM borate buffer (pH10.2)		
	WashSpeed = $25.000 [\mu L/s]$	ReagentBVial: OPA reagent (0.8 mL 400 mM borate buffer + 0.2 mL 75 mg/L OPA + 2 µL 3-mercaptopropionic acid) ReagentCVial: 2.5 mg/mL FMOC ReagentDVial: 50 mM phosphate buffer (pH 3.0)		
	LoopWashFactor = 2.000			
	PunctureOffset = 0.0 [mm]			
	InjectMode = UserProg			
	ReagentAVial = RA1			
	ReagentBVial = RB1			
	ReagentCVial = RC1			
	ReagentDVial = RD1			
	SyringeSpeed = GlobalSpeed			
	SampleHeight = GlobalHeight			
1	UdpDraw From = ReagentAVial, Volume = 50.000	Sample vial contains air sample extractions or standard solutions. Derivatize primary amines using OPA reagent.		
2	UdpMixWait Duration = 1			
	UdpDraw From = ReagentBVial,			
3	Volume = 2.500			
4	UdpMixWait Duration = 1			
5	UdpMixNeedleWashVolume = 100.000			
6	UdpDraw From=ReagentAVial, Volume=2.500			
7	UdpDraw From = SampleVial, Volume = 1.000, SampleHeight = 5.000			
8	UdpDraw From = Air, Volume = 6.000			
9	UdpMoveSyringe Load = 6.000 UdpMoveSyringe Unload = 6.000	Repeat Step 9 five times to complete the derivatization.		
10	UdpMixWait Duration = 60			
11	UdpMixNeedleWash Volume = 100.000			
12	UdpDraw From = ReagentCVial, Volume = 1.000,	Derivatize primary and secondary amines using FMOC reagent.		
13	UdpMixWait Duration=1			
14	UdpDraw From = Air, Volume = 7.000			
15	UdpMoveSyringe Load = 7.000 UdpMoveSyringe Unload = 7.000	Repeat Step 15 eight times to complete the derivatization.		
16	UdpMixNeedleWash Volume=100.000			
17	UdpDraw From = ReagentDVial, Volume = 14.000	Adjust the pH value of the solution in the sample loop to match the		
18	UdpMixWait Duration = 1	mobile phase requirement.		
	UdpDraw From = Air, Volume = 15.000			
20	UdpMoveSyringe Load = 15.000 UdpMoveSyringe Unload = 15.000	Repeat Step 20 five times to thoroughly mix the solution.		
21	UdpInjectValve Position = inject			

Results and Discussion

To ensure that the derivatization of primary and secondary amines taking place in the sample loop was completed, the amounts of OPA and FMOC reagents were investigated. Experiments showed that the responses of 2 µg/mL amine standards using OPA (2.5 µL) and FMOC (1 µL) reagents were the same as those using 1/50 amount of OPA and FMOC reagents. This result demonstrated that the amounts of OPA (2.5 µL) and FMOC (1 µL) exceeded the amount required for derivatization of 2 µg/mL amine standards and were enough for the derivatization of common air samples with amine content up to the µg/mL level.

Similar to using a basic buffer mobile phase to separate OPA- and FMOC-derivatized amino acids by reversed-phase HPLC,²⁵ the use of a basic buffer mobile phase for the separation of derivatized aliphatic amines may also produce higher resolution, lower detection limits, and better method reproducibility. Therefore, the Acclaim™ PA2 column that provides enhanced hydrolytic stability from pH 1.5–10 was chosen for use with a basic buffer mobile phase. Figure 2 illustrates good separation of seven amine standards automatically derivatized using OPA (for ethanolamine, methylamine, ethylamine, *n*-propylamine, *n*-butylamine, and *n*-hexylamine) and FMOC (for diethylamine).

Method precision was estimated by making five consecutive injections of a calibration standard with a concentration of 1 mg/L for each amine. The retention time and peak area reproducibilities are summarized in Table 2. Calibration linearity for fluorescence detection of ethanolamine, methylamine, ethylamine, diethylamine, *n*-propylamine, *n*-butylamine, and *n*-hexylamine was investigated by making three consecutive injections of a mixed standard prepared at five different concentrations (i.e., 15 total injections).

The external standard method was used to establish the calibration curve and quantify the analytes in the air samples. Excellent linearity was observed from 0.05 to 5 mg/L when plotting the concentration versus the peak area, and the correlation coefficients were \geq 0.97 for all. The method detection limits (MDL) of each analyte for fluorescence detection, calculated using signal-to-noise = 3, were all \leq 20 µg/L.

Figure 3 shows the chromatograms of an air sample as well as a mixture of ethanolamine, methylamine, ethylamine, diethylamine, *n*-propylamine, *n*-butylamine, and *n*-hexylamine standards. No detectable-level aliphatic amines were found except for diethylamine (Peak 4). The calculated concentration of the peak labeled as diethylamine is 20 µg/L, which is similar to the estimated MDL of diethylamine, indicating the presence of diethylamine in the air sample. A small peak with retention time near that of *n*-propylamine (Peak 5) was found in the air sample. However, comparison of the retention times reveals that this peak was from the blank (water). Recoveries for each aliphatic amine standard in the sample ranged from 85 to 95%, demonstrating good accuracy of the HPLC method.

Column: Acclaim PA2, 3 µm (3.0 × 150 mm)
Mobile Phase: A: 50 mM phosphate buffer, pH 8.5

B: CH₃CN/CH₃OH (1:1, v/v)

Gradient: B: 0–4 min, 35%; 4–7 min, 35–55%; 7–15 min, 55% Flow Rate: 0.6 mL/min

Inj. Volume: 2 µL
Temperature: 40 °C
Detection: Fluores

Λ

-50,000,000

Fluorescence Peaks: 1. Ethanolamine Ex. 230 nm, Em. 450 nm 2. Methylamine 3. Ethylamine

8

Minutes

12

15

500,000,000 7 3 4. Diethylamine 5. *n*-Propylamine 6. *n*-Butylamine 7. *n*-Hexylamine 8. FMOC reagent 5

Figure 2. Chromatograms of aliphatic amine standards (1 μ g/mL each) automatically derivatized using OPA and FMOC.

Table 2. Reproducibility for peak retention time and area

Analyte	Retention Time RSD	Peak Area (RSD)
Ethanolamine	0.16	2.02
Methylamine	0.26	4.16
Ethylamine	0.44	3.04
Diethylamine	0.38	3.23
<i>n</i> -Propylamine	0.49	0.87
n-Butylamine	0.24	2.27
n-Hexylamine	0.08	0.28

Column: Acclaim PA2, $3 \mu m (3.0 \times 150 mm)$ Mobile Phase: A: 50 mM phosphate buffer, pH 8.5

B: CH₂CN/CH₂OH (1:1, v/v)

Gradient: B: 0-4 min, 35%; 4-7 min, 35-55%; 7-15 min, 55%

| Flow Rate: 0.6 mL/min | Inj. Volume: 2 µL | Temperature: 40 °C | Fluorescence | Ex. 230 nm, Em. 450 nm

C-grams: a. Mixture of amine standards c. Air sample

(1 µg/mL each) d. Air sample spiked with a mixture of amine standards (1 µg/mL each)

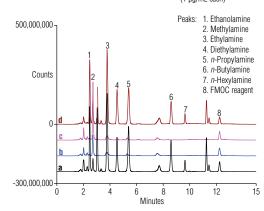


Figure 3. Chromatograms of an air sample and the same sample spiked with a mixture of amine standards.

Conclusion

The work shown here describes an efficient HPLC method using fluorescence detection coupled with automatic precolumn derivatization for the determination of ethanolamine, methylamine, ethylamine, diethylamine, *n*-propylamine, *n*-butylamine, and *n*-hexylamine in air samples. The application of automatic precolumn derivatization reduces the labor required for the determination of aliphatic amines by reversed-phase HPLC

References

- 1. Namiesnik, J.; Jastrzebska, A.; Zygmunt, B. Determination of Volatile Aliphatic Amines in Air by Solid-Phase Microextraction Coupled with Gas Chromatography with Flame Ionization Detection. *J. Chromatogr.*, A 2003, 1016, 1–9.
- Santagati, N.A.; Bousquet, E.; Spadaro, A.; Ronsisvalle, G. Analysis of Aliphatic Amines in Air Samples by HPLC with Electrochemical Detection. *J. Pharm. Biomed. Anal.* 2002, 29, 1105–1111.
- 3. Akyuz, M. Simultaneous Determination of Aliphatic and Aromatic Amines in Indoor and Outdoor Air Samples by Gas Chromatography–Mass Spectrometry. *Talanta* 2007, 71, 486–492.
- Akyuz, M. Simultaneous Determination of Aliphatic and Aromatic Amines in Ambient Air and Airborne Particulate Matters by Gas Chromatography–Mass Spectrometry. *Atmospheric Environment* 2008, 42, 3809–3819.
- Akyuz, M.; Ata, S. Simultaneous Determination of Aliphatic and Aromatic Amines in Water and Sediment Samples by Ion-Pair Extraction and Gas Chromatography–Mass Spectrometry. *J. Chromatogr.*, A 2006, 1129, 88–94.
- 6. Mishra, S.; Singh, V.; Jain, A.; Verma, K.K. Simultaneous Determination of Ammonia, Aliphatic Amines, Aromatic Amines and Phenols at μg/L Levels in Environmental Waters by Solid-Phase Extraction of Their Benzoyl Derivatives and Gas Chromatography— Mass Spectrometry. Analyst (Cambridge, U.K.) 2001, 126, 1663–1668.
- 7. Zhao, X.E.; Li, Y.L.; You, J.M.; Liu, Y.J.; Suo, Y.R. Precolumn Derivatization–High-Performance Liquid Chromatography for the Determination of Aliphatic Amines with Fluorescence Detection and Mass Spectrometric Identification. *Chin. J. Anal. Chem.* 2007, 35, 779–785.
- You, J.M.; Song, C.H.; Yan, T.; Sun, Z.W.; Li, Y.L.; Suo, Y.R. An Improved Reagent for Determination of Aliphatic Amines with Fluorescence and Online Atmospheric Chemical Ionization–Mass Spectrometry Identification. *Anal. Chim. Acta* 2010, 65, 898–905.
- 9. Gao, C.X.; Krull, I.S. Determination of Aliphatic Amines in Air by OnLine Solid-Phase Derivatization with HPLC–UV/FL, *J. Chromatogr. Sci.* **1990**, 28 (3), 102–108.

- Zhang, N.; Zhao, Y.Z.; Zhang, H.S.; Wang, H. Sensitive Determination of Aliphatic Amines by High-Performance Liquid Chromatography with a New Fluorogenic Probe 3-(4-Fluorinebenzoyl)-2-Quinoline Carboxaldehyde. J. Sep. Sci. 2008, 31, 38–46.
- Liu, X.; Wang, H.; Liang, S.C.; Zhang, H.S.
 Determination of Primary and Secondary Aliphatic
 Amines by N-hydroxysuccinimidyl-4,3,2 Naphthapyrone-4-Acetate and Reversed Phase
 High-Performance Liquid Chromatography. *Anal. Chim. Acta* 2001, 441, 45–52.
- Lamba, S.; Pandit, A.; Sanghi, S.K.; Gowri, V.S.; Tiwari, A.; Baderia, V.K.; Singh, D.K.; Nigam, P. Determination of Aliphatic Amines by High-Performance Liquid Chromatography–Amperometric Detection after Derivatization with Naphthalene-2,3-Dicarboxaldehyde. *Anal. Chim. Acta* 2008, 614, 190–195.
- Santagati, N.A.; Bousquet, E.; Spadaro, A.;
 Ronsisvalle, G. Analysis of Aliphatic Amines in Air
 Samples by HPLC with Electrochemical Detection. J.
 Pharm. *Biomed. Anal.* 2002, 29, 1105–1111.
- Hermans, C.; Jonkers, A.C.A.; Bokx, P.K.
 Determination of Amines in the Presence of Excess Ammonia by Ion Chromatography–Mass Spectrometry. J. Chromatogr. Sci. 2010, 48, 544–548.
- Mori, M.; Murad, I.H.H.; Xu, Q.; Hu, W.Z.; Ikedo, I.; Ding, M.Y.; Taoda, H.; Tanaka, K. Vacancy Ion-Exclusion/Adsorption Chromatography of Aliphatic Amines on a Polymethacrylate-Based Weakly Basic Anion-Exchange Column.

 Chromatogr., A 2004, 1039, 129–133.
- Casella, T.G.; Palladino, G.A.; Contursi, M.
 Determination of Aliphatic Amines by Cation Exchange Chromatography with Suppressed
 Conductivity Detection After Solid Phase Extraction.
 J. Sep. Sci. 2008, 31, 3718–3726.
- 17. Hajós, P.; Horváth, K.; Conca, R.; Sarzanini, C. Histidine as a Dipolar Eluent in Ion Chromatography of Aliphatic Amines. *Chromatographia* **2002**, *56*, 103–107.
- 18. Chang, I.H.; Lee, C.G.; Lee, D.S. Development of an Automated Method for Simultaneous Determination of Low Molecular Weight Aliphatic Amines and Ammonia in Ambient Air by Diffusion Scrubber Coupled to Ion Chromatography. *Anal. Chem.* 2003, 75, 6141–6146.
- IonPac CS14 Cation-Exchange Column; LPN 034890-01; Thermo Scientific (formerly Dionex): Sunnyvale, CA, 1997; Documents, Data Sheets; Columns and Accessories Data Sheets; [Online] www.dionex.com/en-us/webdocs/4233-Cs14.pdf (accessed April 29, 2012).

- 20. Application Note 182: Determination of Biogenic Amines in Alcoholic Beverages by Ion Chromatography with Suppressed Conductivity and Integrated Pulsed Amperometric Detections. 2007. Thermo Scientific (formerly Dionex): Documents, Applications. [Online] www.dionex.com/en-us/webdocs/56196-AN182_IC_BiogenicAmines_Alcohol_14Aug2007_LPN1888_02.pdf (accessed April 29, 2012).
- 21. Application Note 183: Determination of Biogenic Amines in Fermented and Non-Fermented Foods Using Ion Chromatography with Suppressed Conductivity and Integrated Pulsed Amperometric Detections. 2007. Thermo Scientific (formerly Dionex): Documents, Applications. [Online] www. dionex.com/en-us/webdocs/57922-AN183_IC_BiogenicAmines_Food_23Aug07_LPN1889_02.pdf (accessed Mar. 27, 2012).
- 22. Saeai, T.; Kato, M.; Teshima, N. Simultaneous Spectrophotometer Analysis of Aliphatic Amines Utilizing Thermochromism of Charge-Transfer Complexes with Tetrabromophenolphthalein Ethyl Ester. *Anal. Sci.* 2005, 21, 1557–1560.
- 23. Molin-Martínez, Y.; Campins-Falco, P.; Herraez-Hernadez, R. A Method for the Determination of Dimethylamine in Air by Collection on Solid Support Sorbent with Subsequent Derivatization and Spectrophotometric Analysis. *J. Chromatogr.*, A 2004, 1059, 17–24.

- 24. Fekete, A.; Malik, A.K.; Kumar, A.; Schmitt-Kopplin, P. Amines in the Environment. *Anal. Chem.* **2010**, 40, 102–121.
- 25. Deng, W.H.; Shen, Y.B.; Li, Z.Y.; Zhang, S.J.; Liu, X.P. Determination of Gamma-Aminobutyric Acid and Seventeen Free Amino Acids in Pinus Massoniana Needles by Reversed-Phase High-Performance Liquid Chromatography with Precolumn Derivatization. *Acta Eco. Sinica* 2008, 28, 4023–4029.
- Hong Yang, H.; Yu, J.Z.; Ho, S.S.H.; Xu, J.H.; Wu, W.S.; Wan, C.H.; Wang, X.D.; Wang, X.R.; Wang, L.S. The Chemical Composition of Inorganic and Carbonaceous Materials in PM2.5 in Nanjing, China. *Atmospheric Environment* 2005, 39, 3735–3749.

www.thermofisher.com/dionex

©2016 Thermo Fisher Scientific Inc. All rights reserved. Sigma-Aldrich is a registered trademark of Sigma-Aldrich Co. LLC. Millex is a trademark of Millipore Corporation.All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details. This information is presented as an example of the capabilities of Thermo Fisher products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others.

Australia +61 3 9757 4486 Austria +43 1 616 51 25 Benelux +31 20 683 9768 +32 3 353 42 94 Brazil +55 11 3731 5140 China +852 2428 3282

Denmark +45 36 36 90 90

France +33 1 39 30 01 10

Germany +49 6126 991 0

India +91 22 2764 2735

Ireland +353 1 644 0064 Italy +39 02 51 62 1267 Japan +81 6 6885 1213 Korea +82 2 3420 8600 Singapore +65 6289 1190 Sweden +46 8 473 3380 Switzerland +41 62 205 9966 Taiwan +886 2 8751 6655 UK +44 1276 691722 USA and Canada +847 295 7500



Part of Thermo Fisher Scientific