

高分解能 GC/Q-TOF による 消防士用防護製品の燃焼生成物の分析



著者

Christiane Hoppe-Jones¹, Shawn Beitel¹, Sofia Nieto², Nathan Eno², Craig Marvin³, and Shane Snyder¹

- ¹ Department of Chemical and Environmental Engineering, University of Arizona, Tucson, AZ,
- ² Agilent Technologies, Inc. Santa Clara, CA
- ³ Agilent Technologies, Inc Wilmington, DE

概要

癌は、消防士の疾病率や死亡率の主な要因の1つです[1,2]。消防士は消火中、煙に加えて、家具、 カーペットなどから発生する、難燃剤などの高い濃度の有機化学物質やその燃焼生成物などにさら されます。フルスペクトル採取モードの高分解能 MS は、複雑なマトリックス中の環境汚染物質のノ ンターゲット分析、特に微量化合物の同定に対して非常に有効です。多環芳香族炭化水素 (PAH)、ポ リ臭化ジフェニルエーテル (PBDE)、その他の燃焼生成物を同定するために、高分解能 Agilent 7250 GC/0-TOF を使用しました。





実験方法

サンプル収集と抽出メソッド



図 1. サンプル収集と抽出ワークフロー

分析

GC/MS 分析には、Agilent 7890B GC システム、新しい高分解能 (m/z 272 に おいて 25,000) Agilent 7250 GC/Q-TOF、および Low-Energy 対応 EI イオン源 を組み合わせて使用しました。

表1に、分析条件を示します。

表 1. Agilent 7250 GC/Q-TOF 分析条件

パラメータ	設定
カラム	Agilent DB-5MS∖ 0.25 mm × 30 m∖ 0.25 µm
注入量	1 µL
注入モード	スプリットレス
スプリット/スプリットレス注入口温度	280 ° C
オーブン温度プログラム	50°Cで3分、 10°C/minで300°Cまで上昇、 7分維持
キャリアガス	ヘリウム、1.5 mL/min、定流量
トランスファライン温度	300 ° C
イオン化モード	標準 El、70 eV Low- Energy El、15 eV および 12 eV
イオン源温度、 70 eV/15 eV 以下	240 ° C/200 ° C
四重極温度	150 ° C
質量範囲	$50\sim$ 1,200 m/z
スペクトル採取レート	5 Hz



図 2. PBDE の分解、およびダイオキシンとフラン構造の形成

データ解析

データの取り込みには、Agilent 7250 精密質量高分解能 GC/0-TOF シス テムを使用しました。最初に、Unknowns Analysis B.08.00 の SureMass 検 出アルゴリズム機能を用いてデータを処理しました。最初の化合物同定 は NIST14 EI ライブラリとのスペクトル比較により実施し、可能な場合は リテンションインデックス (RI) 照合により確認しました (図 3)。Unknowns Analysis の結果を基に、Agilent MassHunter Quantitative Analysis ソフトウェ ア B.08.00 で Quant メソッドを作成し、PAH および PAH 様化合物を半定 量しました (図 4)。未知臭素化化合物の PBDE および分子イオンを、Low-Energy El 化合物スペクトルにより同定しました。MassHunter Qualitative Analysis B.08.00 の Molecular Formula Generator (MFG) を用いて、*m/z* の全同 位体クラスタ、相対同位体アバンダンス、および同位体比を評価し、分子 イオンを確認しました。



図 3. Unknowns Analysis B.08.00 および NIST14 とのスペクトル比較

	17 (B) (M)	Ne (1996, South)		Compo	2N	Mass Accuracy	M7	A	
Sargin						mass Accordcy	191dai	111	
Name FF Nack 365 pror-1	Eals File FF Neck 301 prov-1.D	Type La	VET 2016 15 31 AM	RT 25.207	3.35	-1.8980	216.0930	1	
FF_beck_X8_peop-J FF_beck_X8_peop-3	FF Neck NB pror 3.0 FF Neck NB pror 3.0	Sanple Sanple	12/12/16/11 11 AM 12/12/16/11 11 AM	2222	7.01	0.2373	216.0935	1	
FF_Neck_205_point1 FF_Neck_305_point2	FF_Neck_308_post 1.0 FF_Neck_308_post 2.0	Sargle Sargle	1212016 1245 PM	23 229	6.26	-2.4555	216.0929	1	
I Maimat post-1	FF Individ post-10	Sample	157 2016 3 00 PM	222	4.55	-0.1726	216.0934	1	
F plainer, post-3	FF yearer past-3.0	Sarger	10/10016-4 20 PM	27.236	7.73	-0.1259	216.0934	1	
-					7.66	0.2633	216.0935	1	
A I II	IN SE AT	4 6 4	X A A A A	4.6	2.86	-1.2214	216.0931	1	
a pose ta "usuar" he	8011P				0.64	-1.1653	216.0931	1	
					9.70	-1.3967	216.0931	1	
	~~		nm	100					
a nja nja	ฟา ม่น ม	14 22/16 2	1/11 2/22 2/22	23/24	22	ע אלת מלת כלו אלת	and 21.50 21.4 Accuration Tom	25'42 e (Mini	

図 4. Agilent MassHunter Quant B.08 を用いた PAH の定量

結果と考察

PAH の同定

火災後のすべてのサンプル中の PAH を同定しました。火災後のヘルメット上で、50 を超える PAH 種を同定しました。NIST14 を用いて同定した規制対象 PAH は、リテンションインデックスにより確認しました。表2 は、 火災後のヘルメットで同定された PAH および PAH 様化合物の例を示しています。 同定された分子式について可能性のある構造を表 2 に示しています。これらは、火災後の空気中および消防士の皮膚から Fernando 他 [3] によって検出された PAH に類似しています。

表 2. 同定されたPAH を組み合わせたリスト。注: 大部分の PAH および PAH 様分子式において、複数の異性体を同定

化合物/分子式	m/z	平均質量誤差 (ppm)	化合物/分子式	m/z	平均質量誤差 (ppm)	化合物/分子式	m/z	平均質量誤差 (ppm)
規制対象 PAH			その他の PAH			0 含有 PAH		
ナフタレン [C ₁₀ H ₈]	128.0626	1.01	$C_{10}H_{12}$	132.0934	0.91	C ₁₃ H ₈ O	180.057	0.90
アセナフチレン [C ₁₂ H ₈]	152.0626	0.82	$C_{11}H_{14}$	146.109	1.07	$C_{16}H_{10}O$	218.0726	1.47
アセナフテン [C ₁₂ H ₁₀]	154.07825	0.35	$C_{11}H_{10}$	142.0777	0.55	C ₁₇ H ₁₀ O	230.0726	1.99
フルオレン [C ₁₃ H ₁₀]	166.07825	1.15	$C_{12}H_{16}$	160.1247	1.22	C ₁₇ H ₁₀ O ₂	246.0675	1.27
フェナントレン [C ₁₄ H ₁₀]	178.07825	1.25	$C_{12}H_{12}$	156.0934	1.07	C ₁₈ H ₁₀ O	242.0726	2.03
アントラセン [C ₁₄ H ₁₀]	178.07825	0.96	$C_{13}H_{12}$	168.0934	1.14	$C_{18}H_{10}O_{2}$	258.0675	1.15
フルオランテン [C ₁₆ H ₁₀]	202.07825	1.06	$C_{14}H_{14}$	182.109	0.59	C ₂₀ H ₁₂ O	268.0883	1.68
ピレン [C ₁₆ H ₁₀]	202.07825	1.51	$C_{15}H_{12}$	192.0934	0.53	_		
ベンゾ[a]アントラセン [C ₁₈ H ₁₂]	228.0939	1.27	$C_{15}H_{10}$	190.0777	2.57		$\bigcirc \neg _$	
クリセン [C ₁₈ H ₁₂]	228.0939	1.07	C ₁₈ H ₂₂	238.1716	1.07	-	- YA	
ベンゾ[b]フルオランテン [C ₂₀ H ₁₂]	252.0939	1.81	C ₁₈ H ₁₈	234.1403	1.04			
ベンゾ[k]フルオランテン [C ₂₀ H ₁₂]	252.0939	2.23	$C_{17}H_{12}$	216.0934	0.87	U ₁₃ H ₈ U	U ₁₆ H ₁₀ U	
ベンゾ[a]ピレン [C ₂₀ H ₁₂]	252.0939	1.79	$C_{18}H_{10}$	226.0777	0.92		$\wedge \wedge$	1
インデノ[1,2,3 -cd]ピレン [C ₂₂ H ₁₂]	276.0939	1.65	$C_{20}H_{14}$	254.109	1.05	ųμ	YNY	
ジベンズ[a,h]アントラセン [C ₂₂ H ₁₄]	278.10955	1.55	$C_{19}H_{14}$	242.109	0.58	CHO.	C ₁₈ H ₁₀ O	
ベンゾ[ghi]ペリレン [C ₂₂ H ₁₂]	276.0939	1.73	$C_{19}H_{12}$	240.0934	2.17	-17-10-2	10 10	
_			$C_{22}H_{14}$	278.10955	0.91	化合物/分子式	m/z	平均質量誤差 (ppm)
$ \land \land$	(Ω)	$\hat{\Box}$	$C_{24}H_{18}$	306.1403	1.67	N 含有 PAH		
	Ų,	QQ	$C_{22}H_{12}$	276.0939	2.25	C ₁₃ H ₉ N	179.073	1.24





_	$C_{15}H_9N$	203.073
	C ₁₇ H ₁₁ N	229.0886
	C ₁₉ H ₁₁ N	253.0886
		\sim

 $C_{13}H_9N$

 $C_{12}H_9N$



167.073

1.55

1.52

0.63

1.08

相対アバンダンスを基にした分析の結果、火災後のヘルメットで測定され た主な PAH は、フルオランテン、ピレン、および大型の PAH でした。図5 は、火災後のヘルメットの PAH および PAH 様化合物のアバンダンスを示 しています。 図6は、すべてのサンプル中のPAH、0含有PAH、およびN含有PAHの合計を示しています。火災後のヘルメットのPAHは、皮膚での濃度を超えていました。火災後の皮膚サンプルは、すべてのグループのPAHが高い濃度を示していました。



図 5. 火災後のヘルメットの拭き取りで検出された PAH および PAH 様化合物



図 6. 火災前後の拭き取りで同定された総 PAH

臭素化化合物

臭素化化合物は、火災後のサンプルでのみ検出されました。これは、サン プルが煙にさらされたことに起因することを示しています。

火災後の首とヘルメットからは、難燃剤を含む燃焼物質に起因するトリ -BDE、テトラ-BDE、ペンタ-BDE、およびヘキサ-BDEの異性体が検出されま した。表3に相対アバンダンスを示します。リテンションタイムは、未変性 標準を用いて検証しました。 その他の Br 含有化合物は火災後のヘルメットのみから検出され、首から は検出されませんでした。これらの化合物は、PBDE 燃焼生成物 (例えば $C_6H_4Br_20$) および臭素化 PAH (例えば $C_{16}H_9Br$) の可能性があります。図 7 と 図 8 は、異なるイオン化エネルギーでのいくつかの臭素化化合物の質量 スペクトルを示しています。Low-Energy EI は、未知化合物の分子イオン同 定に有用でした。

表 3. 火災後のサンプルで同定された PBDE およびその他の臭素化化合物のリスト

分子式	m/z	質量誤差 (nnm)*	分解能	MFG スコア	標準により 検証	リテンション タイム (分)	首: 火災前、 % アバンダンス	首: 火災後、 % アバンダンス	ヘルメット: 火災後、 % アバンダンス
	2/0 8623	1.7	27226	08.34	17.000	13 15	nd	nd	15.26
G6H4DI2U	249.0023	1.7	27330	90.34		13.15	n.u.	n.u.	10.20
$C_{14}H_{21}BrO$	284.077	0.92	26911	90.92		16.409	n.d.	n.d.	31.17
$C_7H_8Br_2N_2$	277.9049	0.22	27495	88.15		19.891	n.d.	n.d.	5.72
$C_7H_6Br_2N_2O$	291.8841	0.99	26072	90.08		20.832	n.d.	n.d.	13.84
$C_{10}H_9BrN_2$	235.9944	1.27	27136	97.41		21.122	n.d.	n.d.	1.53
$C_{16}H_9Br$	279.9882	1.69	26187	98.71		24.609	n.d.	n.d.	4.39
$C_{15}H_{14}Br_2O_2$	383.9355	0.25	29880	95.36		25.88	n.d.	n.d.	0.32
$C_{12}H_7Br_3O$	403.8047	0.45	31609	90.67	トリ-BDE	22.687	n.d.	1.1	0.29
		0.23	29157	96.13		23.063	n.d.	1.41	0.4
$C_{12}H_6Br_4O$	481.7152	1.36	30425	91.87	テトラ-BDE	24.835	n.d.	0.9	0.4
		2.62	31364	95.40		25.175	n.d.	58.84	13.09
		1.15	27322	93.01		25.507	n.d.	0.69	0.3
$C_{12}H_5Br_5O$	559.6257	2.59	30367	95.05	ペンタ-BDE	26.684	n.d.	8.99	2.74
		2.03	30813	94.57		27.159	n.d.	25.29	9.26
		1.24	29448	91.29		27.974	n.d.	1.1	0.35
$C_{12}H_4Br_6O$	637.5357	0.7	30817	91.38	ヘキサ-BDE	28.335	n.d.	0.83	0.51
		1.7	31299	94.73		29.058	n.d.	0.85	0.43

*質量誤差は、全同位体クラスタの重み付け平均質量誤差として計算



図 7. 火災後の首の拭き取りで検出された PBDE の例。MFG スコアは、質量精度、同位体間隔、および相対同位体アバンダンス比を基にして計算





図 8. 異なるイオン化エネルギーでの未知臭素化化合物スペクトルの例。スペクトルは、MFG ツールと Fragment Formula Annotation を用いてアノテートされ、 アノテートされていないイオンは、共溶出化合物に起因する汚染物質イオンの可能性あり。左側に示したのは、70 eV および Low-Energy EI スペクトルに 基づいた構造候補

結論

火災後に首を拭き取ってみたところ、フードは煙関連の汚染物質を完全 には遮断できていないことがわかりました。火災後の消防士の首とヘル メットから、PAH、PBDE、およびその他の臭素化化合物が検出されました。 臭素化化合物は検出限界が低いため、消防士が煙にさらされたことを示 すバイオマーカーとして機能します。

適用した GC/Q-TOF メソッドは、多数の PAH、PBDE、およびその他の臭素 化燃焼生成物の検出に対して、非常に高感度で選択性があることが証明 されました。

Low-Energy 対応 El イオン源は、未知臭素化化合物の分子イオン同定に有用でした。

参考文献

- M. N. Bates. "Registry-based case-control study of cancer in California firefighters" *Am. J. Ind. Med.* **50** (5), 339-344 (2007).
- D. Kang, et al. "Cancer incidence among male Massachusetts firefighters 1987-2003" Am. J. Ind. Med. 51(5), 329-335 (2008).
- S. Fernando. "Evaluation of firefighter exposure to wood smoke during training exercises at burn houses" *Environ. Sci. Tech.* 50, 1536-1543 (2016).

詳細

本文書のデータは代表的な結果を記載したものです。 アジレント製品とサービスの詳細については、アジレントのウェブサイト www.agilent.com/chem/jp をご覧ください。

ホームページ

www.agilent.com/chem/jp

カストマコンタクトセンタ 0120-477-111 email japan@agilent.com

本製品は一般的な実験用途での使用を想定しており、医薬品医療機器 等法に基づく登録を行っておりません。本文書に記載の情報、説明、 製品仕様等は予告なしに変更されることがあります。

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2017 Printed in Japan, July 6, 2017 5991-8197JAJP

