

下水汚泥における環境汚染物質の ターゲット化合物の定量と サスペクト化合物のスクリーニングの 同時実行

Agilent 6546 LC/Q-TOF

著者

James S. Pyke Agilent Technologies, Inc. Santa Clara CA USA

Gabrielle Black University of California, Davis CA USA

Kai Chen Agilent Technologies, Inc. Santa Clara CA USA

Tarun Anumol Agilent Technologies, Inc. Santa Clara CA USA

Thomas M. Young University of California, Davis CA USA

概要

環境汚染物質のリストは拡大しており、規制はますます厳しくなっています。さまざまな環境機関が、高 速、高精度、高感度の分析ツールを必要としています。環境ラボが取り扱うサンプルには幅広い分析が 必要です。これらは、生態系の保護を目的とした毒性の高い汚染物質の微量分析から、製造時に想定 されていた使用期間を超えて保管されることが多い汚染物質を含む大規模ライブラリのスクリーニング まで、多岐にわたります。データを取り込み、定量して結果をレポートする技法と手順は、通常、ターゲッ ト LC-TQ 技術を用いて実現されてきました。しかし、新しい環境毒性物質は増え続けており、そのリス トをモニタリングすることは、環境科学者にとって依然として課題となっています。

このアプリケーションノートでは、Agilent 6546 LC/Q-TOF MS システムによる定量とスクリーニン グを組み合わせたワークフローを評価します。6546 は、広いダイナミックレンジと高質量分解能を 兼ね備えており、取り込みスピードも優れています。アジレントのパーソナル化合物データベースラ イブラリ (PCDL) の十分に精査された MS/MS スペクトルとリテンションタイム、そして更新された MassHunter Quantitative Analysis ソフトウェアと組み合わせることで、ターゲット化合物のシームレ スな定量と同時に、サスペクト化合物のスクリーニング機能を実行するワークフローが実現します。ソフ トウェアの拡張とハードウェア機能によって、規制化合物の高速かつシンプルな定量が可能になり、一 度の注入で何千もの汚染物質をスクリーニングできます。

はじめに

下水汚泥は、濃縮された化合物の複雑な混合 物で、多くの場合、土壌還元処理されます。土 壌還元された下水汚泥で、一般消費財由来の 残留性有毒化学物質を規制に従ってモニタリ ングするには限界があります。しかし、これら 多数の化合物の毒性値は依然として明らかに されておらず、排出の下流における生態系へ の影響について、リスクを調査し、軽減する必 要があります。課題は、提供される製品の増 加に伴い、毒性物質とその変換生成物のリス トが拡大し続けていることです。一般消費財 由来の残留性化学物質が、堅牢な廃液処理 技術では除去できないとき、これらの化合物 を幅広くスクリーニングすることにより、物質 が下流に与える影響の全体像を明らかにする ことが可能になります。

内分泌かく乱物質は、ホルモンによってコント ロールされる生物系に干渉する化学物質であ るため、下水処理場のプロセスでモニタリン グされています。下水汚泥中の内分泌活性有 機環境汚染物質の分析について説明する分析 メソッド¹を更新し、6546 四重極飛行時間型 (Q-TOF) システムを最大限に活用できるよう にしました。6546 LC/Q-TOF は、広いダイナ ミックレンジと高質量分解能機能を兼ね備え ており、取り込みスピードも優れています。ま た、Agilent 1290 Infinity II 液体クロマトグラ フィー (LC) システムと組み合わせて高速 LC グラジエントを適用し、クロマトグラフィー分 離能を向上させることにより、ハイスループッ ト分析を実施することが可能になります。 これらの機能を組み合わせることで、既知の 毒性物質の高速な定量と、その他多数の存在 する可能性のある毒性物質の存在のモニタリ ングが可能になり、実行した作業の価値が高 まります。さらに、データ非依存型取り込み (DIA)機能により、新しい毒性物質が発見さ れた際には遡及的な解析も実行できます。

今回、選択した化合物を下水汚泥マトリックス にスパイクして、6546 LC/Q-TOF システムの 定量能力を評価しました。下水汚泥中で以前 検出されたことがあり、内分泌物の活性特性 に影響を与えると考えられる化合物約 50 を リストにまとめました。メソッドバリデーション では、これらの化合物のうち広範なリストの 牛理化学的性質を代表する 12 個のサロゲー トを選択しました。スパイクしたサロゲート は、クロマトグラム全体を通して溶出するのに 十分な化学的多様性を備えているため、サロ ゲート化合物のリテンションタイム (RT) を、 同じ化合物を測定する公開済みの分析メソッ ドと相関付けました。次に、モデルを使用して、 同じデータファイルから幅広い範囲の毒性物 質の RT を見積もりました。さらに、RT の相 関性がない化合物もモニタリングしました。

スパイクしたサロゲートに加えてモニタリング された、十分に精査された MS/MS スペクト ルを持つ化合物は合計 4,856 個でした。シン プルなデータ解析ワークフローで、高質量分 解能データに基づく十分に精査された化合物 ライブラリで提供されている、化合物の既知 のプリカーサとフラグメントの質量を抽出しま す。次に、以下の作業を実行します。

- 抽出した質量の質量精度と共溶出の 測定
- 既知の理論的同位体パターンと正確に 測定されたパターンとの比較
- 既知または予測 RT と測定値の比較

推定上の同定は、SANTE ガイドライン³の推 奨どおり、基本的な同定基準に従っています。 また、ソフトウェアでは確認プロセスに重点を 置き、偽陽性の可能性を低減しています。従 来の定量プロセス (LC/TQ ワークフローと類 似) でスパイクしたサロゲートを検出して定量 することでワークフローを確認した後、十分 に精査された化合物データベースで新しい毒 性物質をさらに拡張することにより、存在する 可能性のある汚染物質を遡及的に検出してい ます。

実験方法

下水汚泥サンプルのサンプル前処理手順を故 意にできるだけ非化学的に選択することによ り、可能な限り幅広い範囲の環境汚染物質を 検出できます。高品質の溶媒と消耗品を使用 して、誤った結果を引き起こす可能性のある 化合物が導入されないようにしました。

試薬および調製

溶媒は、Agilent LC/MS グレードのアセトニ トリル (p/n G2453-85050)、 ギ酸 (Merck、 5330020050)、およびフッ化アンモニウム (Sigma-Aldrich、338869) により調製しま した。リファレンスマス溶液は、100 µL の Agilent HP-0921 および 200 µL のプリン (p/n G1969-85001) を 100 mL の 5 % 水 アセトニトリル溶液に加えて前処理しました。 ESI-L キャリブラント溶液は、機器のマニュア ルに従って調製しました。標準は、カリフォル ニア大学デービス校において 20 % メタノール 水溶液で前処理しました。

標準とサンプル前処理

表1に、ターゲット化合物のリストを示します。 6546 Q-TOF LC/MS システムのダイナミック レンジと感度を評価するために、標準をキャリ ブレーションレベル 1,000、750、500、250、 100、50、25、10、5、2.5、1、0.5、0.25、お よび 0.1 ppb において 20 % メタノール水溶 液で調製しました。データを正規化するため の内部標準は使用しませんでした。サンプル は米国カリフォルニア州にある水処理施設か ら入手して、既報に従って前処理し¹、ターゲッ ト化合物をサンプル前処理前後に各々 200 ppb でスパイクして、回収率とサンプルのマト リックス抑制効果を評価しました。

装置構成

表 2 に示すモジュールで構成された Agilent 1290 Infinity II LC システムを、6546 Q-TOF LC/MS システム (G6546A) と組み合 わせました。

表1. ターゲット化合物

名前	CAS	分子式	中性質量
AHTN/トナリド (フィクソリド)	1506-02-01	C ₁₈ H ₂₆ O	258.19837
カルバマゼピン	298-46-4	C ₁₅ H ₁₂ N ₂ O	236.09496
DEET/ジエチルトルアミド	134-62-3	C ₁₂ H ₁₇ NO	191.13101
ジクロフェナク	15307-86-5	C ₁₄ H ₁₁ Cl ₂ NO ₂	295.01668
ジヒドロジャスモン酸、メチルエステル	24851-98-7	C ₁₃ H ₂₂ O ₃	226.15690
エファビレンツ	154635-17-3	$C_{14}H_9CIF_3NO_2$	315.02739
フルニキシン	38677-85-9	$C_{14}H_{11}F_3N_2O_2$	296.07726
フルオキセチン	54910-89-3	C ₁₇ H ₁₈ F ₃ NO	309.13405
フルボキサミン	54739-18-3	$C_{15}H_{21}F_3N_2O_2$	318.15551
ラモトリギン	84057-84-1	C ₉ H ₇ Cl ₂ N ₅	255.00785
メフェナム酸	61-68-7	C ₁₅ H ₁₅ NO ₂	241.11028
メトプロロール	37350-58-6	C ₁₅ H ₂₅ NO ₃	267.18344
ミコナゾール	22916-47-8	C ₁₈ H ₁₄ Cl ₄ N ₂ O	413.98602
ノルゲストレル	797-63-7	C ₂₁ H ₂₈ O ₂	312.20893
スルファメトキサゾール	723-46-6	C ₁₀ H ₁₁ N ₃ O ₃ S	253.05211
トリクロカルバン	101-20-2	C ₁₃ H ₉ Cl ₃ N ₂ O	313.97805
トリメトプリム	738-70-5	C ₁₄ H ₁₈ N ₄ O ₃	290.13789
エストロン (E1)	53-16-7	C ₁₈ H ₂₂ O ₂	270.16198
エチニルエストラジオール (EE2)	57-63-6	C ₂₀ H ₂₄ O ₂	296.17763
2-フェニルフェノール (オルトフェニルフェノール)	90-43-7	C ₁₂ H ₁₀ O	170.07316
ゲムフィブロジル	25812-30-0	C ₁₅ H ₂₂ O ₃	250.15689
エストリオール	50-27-1	C ₁₈ H ₂₄ O ₃	288.17254
4-tert-オクチルフェノール (4-(1,1,3,3-テトラメチルブチル)フェノール)	140-66-9	C ₁₄ H ₂₂ O	206.16707

表 2. Agilent 1290 Infinity II UHPLC の条件

モジュール	ポジティブイオン化パラメータ	ネガティブイオン化パラメータ			
	溶媒 A1) 0.1 % ギ酸水溶液 溶媒 B1) 0.1 % ギ酸アセトニトリル溶液	溶媒 A2) 1 mM フッ化アンモニウム水溶液 溶媒 B2) アセトニトリル			
	流量: 0.4 mL/min 最大圧力上限: 1,300 bar (動作圧力は 450 bar 未満)				
ハイスピードポンプ (G7120A)	グラジエント: 時間(分) %B 0.00 2.00 0.50 2.00 15.50 100.00 19.50 100.00 20.00 2.00 ストップタイム: 20.00 分 ポストタイム: 1.00 分				
マルチサンプラ (G7167B)	注入量: 1 µL マルチウォッシュ:シートバックフラッシュおよびニードル洗浄、100 % イソプロパノール、 100 % アセトニトリル、100 % 水それぞれを 5 秒				
マルチカラムサーモスタット (G7116B)	カラム温度: 30°C カラム: Agilent InfinityLab Poroshell EC-C18、2.1 × 100 mm、1.9 µm (p/n 685775-924)				

表 2 に示したパラメータにより、図 1A に表さ れる LC グラジエントが生成されました。下水 汚泥、ブランク、およびキャリブレータの 160 回注入によるポンプ圧力曲線を重ねて表示し ています (図 1B)。ポジティブイオン化モード (赤) とネガティブイオン化モード (黒) での代 表的なトータルイオン (TIC) クロマトグラムも 示します (図 1C)。

表 3 は、6546 LC/Q-TOF システムにおいて DIA モードで化合物を測定するのに使用した 条件を示しています。



図 1.1290 Infinity II LC システムに注入された下水汚泥のクロマトグラフィー結果

表 3. 6546 LC/Q-TOF LC/MS システム (G6546A) の条件

パラメータ	ポジティブイオン化パラメータ	ネガティブイオン化パラメータ					
Agilent Jet-Stream イオン源							
ドライガス温度	225						
ドライガス流量	12						
ネブライザ	30						
シースガス温度	350						
シースガス流量	350						
キャピラリー電圧	3500						
ノズル電圧	500						
フラグメンタ電圧	110						
	チューンモード						
イオン極性	ポジティブ	ネガティブ					
質量範囲	低 (1,700 m/z)						
スライサモード	高分解能						
	取り込みモード						
	50 ~ 1,050 m/z						
スピード	8 スペクトル/秒						
コリジョンエネルギー	0、10、20、40 V						
参照質量補正	ボトルA使用を有効化						
参照質量	121.050873 (M+H)* ブリン付加物 119.03632 (M-H) ⁻ ブリン付加物 922.009798 (M+H)* HP-0921 付加物 940.001473 (M+F) ⁻ HP-0921 付						

図 2 は、ターゲット化合物の抽出イオンクロマ トグラム (EIC) を示しています。

データ解析ワークフロー

自動化ルーチンワークフローにおいて Agilent MassHunter Acquisition (バージョ ン10.0.111) を使用し、ポジティブおよびネガ ティブイオン化モードで All Ions MS/MS デー タを取り込みます。データは MassHunter Quantitative Analysis (TOF 用) (バージョン 10.1 プロトタイプ) で Agilent SureMass 技 術³により自動的に処理されるため、ターゲッ ト化合物を高速でより正確に定量し、サスペク ト化合物も同時に検出できます。

MassHunter 定量メソッドは、アジレントの十 分に精査されたパーソナル化合物データベー スライブラリ (PCDL) から化合物をインポート して設定します。Agilent PCDL でのデータの 品質確認と、将来出現する新規汚染物質を追 加する推奨プロセスについて、概要が示され ています⁴。プリカーサイオンに対して定量イ オンを設定し、各化合物に対して最低 2 個の MS/MS フラグメントイオンを確認イオンとし て設定しました。

アジレントの十分に精査された環境水スク リーニング PCDL には、以前に説明した分析 メソッド⁵の RT を厳選して記載しています。図 4 のベン図に内容をまとめて示します。



図2.表1で定義されたターゲット化合物の抽出イオンクロマトグラム





図3.ターゲット化合物の定量とサスペクト化合物のスクリーニングのワークフロー



図4.水スクリーニングのパーソナル化合物データベースライブラリの内容

このアプリケーションで使用されている LC メ ソッドは、環境水スクリーニング PCDL で RT を精査するために使用されているメソッドとは 異なるため、両方の分析で共通のターゲット 化合物を使用して、この分析でのサスペクト化 合物の RT をモデル化しました。図 5 は、両 方のメソッドで RT が既知の 9 個の化合物 (黒 の円)、原点に適合させた指数曲線、および このアプリケーションノートで説明されている LC メソッドを用いて分析した際の、環境水ス クリーニング PCDL の化合物の予測 RT を示 しています。予測 RT (青の円)を、サスペクト 化合物を検索した RT ウィンドウを表すエラー バーとともに示しています。

結果と考察

ターゲット化合物の定量能力

下水汚泥中にスパイクしたサロゲートをモニ タリングすることにより、6546 LC/Q-TOF シ ステムに適用した分析メソッドの定量能力を 評価しました。表 4 に示した化合物の直線ダ イナミックレンジは、0.1~1,000 ppb の範 囲のキャリブレーション標準の直線回帰によ り評価しました。S/N比が3未満のキャリブ レーションレベルは除去しており、検出下限 (LLOD) を表 4 にレポートしています。 最低の キャリブレーションレベルでは S/N ≥ 3 でし た。表4に示した飽和限界(LOS)は直線ダイ ナミックレンジの上限です。高濃度のキャリブ レーションポイントは、決定係数 (R²) が 0.99 を超えるまで除去し、直線を 1/x の重み付け (ここで、x は濃度) で適合させたときの各キャ リブレータの定量精度は ±20% 未満でした。



図 5. 分析メソッド間で共通のサロゲートのリテンションタイムの相関

表 4.6546 LC/Q-TOF の定量能力

			LLOD	LOS
名前	定量イオン	R ²	(ppb)	(ppb)
AHTN/トナリド (フィクソリド)	(M+H)*	> 0.99	5.0	1,000
カルバマゼピン	(M+H)*	> 0.99	1.0	250
DEET/ジエチルトルアミド	(M+H)*	> 0.99	0.50	500
ジクロフェナク	(M+H)*	> 0.99	25	> 1,000
ジヒドロジャスモン酸、メチルエステル	(M+H)+	> 0.99	50	> 1,000
エファビレンツ	(M+H)*	> 0.99	5.0	> 1,000
フルニキシン	(M+H)*	> 0.99	0.50	2,500
フルオキセチン	(M+H)*	> 0.99	0.50	> 1,000
フルボキサミン	(M+H)*	> 0.99	1.0	> 1,000
ラモトリギン	(M+H)*	> 0.99	0.10	100
メフェナム酸	(M+H)*	> 0.99	5.0	> 1,000
メトプロロール	(M+H)*	> 0.99	0.50	> 1,000
ミコナゾール	(M+H)*	> 0.99	0.50	500
ノルゲストレル	(M+H)*	> 0.99	2.5	750
スルファメトキサゾール	(M+H)*	> 0.99	50	> 1,000
トリクロカルバン	(M+H)*	> 0.99	50	> 1,000
トリメトプリム	(M+H)*	> 0.99	0.10	100
エストロン(E1)	(M-H) ⁻	> 0.99	2.5	> 1,000
エチニルエストラジオール (EE2)	(M-H) ⁻	> 0.99	5.0	> 1,000
2-フェニルフェノール (オルトフェニルフェノール)	(M-H) ⁻	> 0.99	25	> 1,000
ゲムフィブロジル	(M-H) ⁻	> 0.99	5.0	> 1,000
エストリオール	(M-H) ⁻	> 0.99	5.0	> 1,000
4-tert-オクチルフェノール (4-(1,1,3,3-テトラメチルブチル)フェノール)	(M-H) ⁻	> 0.99	5.0	> 1,000

大半のケースで、線形適合に 1,000 ppb の キャリブレーションサンプルを含んでも LOS は観察されませんでした。ただし、化合物を 重み付け 1/x (ここで、x は濃度)、非線形指 数曲線回帰で適合させた際、ラモトリギン(効率的にイオン化されたと思われる化合物)以外のすべての場合において、より高いキャリブレーションポイントが含まれており、0.5 μLの

注入時に化合物を最大 1,000 ppb で正確に 定量できました。図 6 は、この分析でのター ゲット化合物の検量線を示しています。



図 6. ターゲット化合物の検量線



図 6. ターゲット化合物の検量線 (続き)

ラモトリギンの場合、検出下限 (LLOD) は観 察されませんでした。これは、最低のキャリブ レータ (0.1 ppb) を注入した際、S/N > 3 の ピークが観察されたためです。この分析メソッ ドでは 1.0 μ L の注入量を使用したため、濃 度 100 ppb はカラムの注入では 100 pg と 同等です。 ターゲット化合物の結果を、LC-TQ データ と同じ方法で示します (図 7 を参照)。クオン ティファイアの積分と予想される RT (図 7A) および確認イオンの共溶出 (図 7B、キャリブ レータから決定された予想比に応じてスケー リング) は、LC/TQ および LC/Q-TOF 取り 込みメソッド間で共通です。精密質量測定で の小数点以下の測定値および予想同位体パ ターン (図 7C、赤のボックス) と、測定した 同位体パターン(図7C、黒のスペクトル)を 比較する機能により、既知の化学式と天然の 同位体アバンダンスが指定された場合、化合 物同定において非常に高いレベルの信頼性 が達成されます。



図 7.500 ppb でスパイクした下水汚泥中のラモトリギンの結果

サスペクト化合物のスクリーニング

LC Screener ツールを、SANTE ガイドライン ²の基準に従って MassHunter Quantitative ソフトウェアの推定同定機能カラーコードに 組み込みました。図 8 の緑色は、2 つを超え るイオン (プリカーサイオンやフラグメントイオ ン) が必要な質量精度で測定され、共溶出し、 予想された RT 範囲 (既知の場合) に収まって いることを示しています。さらに、プリカーサ イオンの同位体パターンも確認しました。図 8 に示すように、ネガティブイオン化モードで測 定されることが想定された 6 個のターゲット 化合物をすべて確認しました。2 個の追加の 化合物もネガティブイオン化モードで確認しま した。オレンジは確認が必要な化合物を示し ており、赤は選択したサンプルで化合物が検 出されなかったことを示しています。



図8. サスペクト化合物の推定上の同定に使用した基準を示す LC Screener ツール

下水汚泥中にスパイクしたサロゲートを検出 して、データ解析ワークフローを確認した後、 スパイクした化合物と他の推定同定したネガ ティブイオン化化合物をレポートしています。 図 9 は化合物を抜粋したものを示していま す。 ネガティブおよびポジティブイオン化化合 物の両方に関する詳細なレポートを表 4 にま とめて示します。 表 4. スパイクした下水汚泥サンプル中で確認済み、 確認が必要、および検出されなかった化合物の まとめ

イオン化 モード	確認済み	確認が 必要	検出 されず
ポジティブ	18	159	3,998
ネガティブ	8	47	1,267

		Scree	ining S	ing Summary Report				Aglient Trusted Answers		
Sample	name: NEG_AllIons	_PstSpk_3		Good	8	Wai	rning 47	Error	1267	
Status	Screening Summary Report	Formula	R.T.	R.T. Diff.	Match Score	Target Ion	Mass Accuracy	# of Qualified Tons	Final Conc.	
1	Xanthohumol	C21H22O5	9.411	9.310		353.1394	-0.61 PPM	3		
1.1	Lauryl hydrogen sulfate	C12H26O4S	8.249	8.144		265.1479	0.11 PPM	2		
1.1	Harmine	C13H12N2O	4.923	4.825		211.0877	-1.28 PPM	3		
1.1	THC / delta9-Tetrahydrocannabinol (Δ 9-THC)	C21H30O2	12.739	12.645		313.2173	-0.57 PPM	3		
1.1	(-)CP-55,940	C24H40O3	8.981	8.890		375.2905	-0.77 PPM	2		
1.1	(±)CP-55,940	C24H40O3	8.981	8.890		375.2905	-0.77 PPM	2		
1	(–)-THC-COOH / (–)-11-Nor-9-Carboxy- tetrahydrocannabinol	C21H28O4	9.488	9.382		343.1915	-0.52 PPM	3		
1.1	4-Methylphenol (p-Cresol)	C7H8O	7.370	7.290		107.0502	7.94 PPM	2		
1.1	Fenofibric acid	C17H15ClO4	5.606	5.498		317.0586	-1.73 PPM	2		
1.1	Bisphenol E	C14H14O2	4.634	4.532		213.0921	0.19 PPM	2		
1	BPS / Bisphenol S	C12H10O4S	4,522	4,415		249.0227	0.02 PPM	4		
	Silibinin	C25H22O10	5.707	5.610		481.1140	-1.58 PPM	3		
	Losartan	C22H23CIN6O	5.482	5.379		421.1549	-0.12 PPM	2		
1 - 1 - 1 - 1 - 1 - 1 - 1 - 1 - 1 - 1 -	Veratramine	C27H39NO2	14,185	14.087		408.2908	-0.96 PPM	2		
į.	BKF (Cyanox 2246) (2,2'-methylene-bis(6-tert- butyl-4-methylphenol))	C23H32O2	12.898	12.796		339.2330	0.43 PPM	2		
1	Curcumin	C21H20O6	8.113	8.011		367.1187	0.45 PPM	2		
1	Iloprost	C22H32O4	6.119	6.037		359,2228	-0.23 PPM	2		
i i	Phenylpyruvic acid	C9H8O3	1.530	1.424		163.0401	1.24 PPM	3		
1	THC-COOH / 11-Nor-9-Carboxy- tetrahydrocannabinol	C21H28O4	9.488	9.382		343.1915	-0.52 PPM	3		
1.1	8-Hvdroxvefavirenz	C14H9CIF3NO3	8.143	8.044		330.0150	0.55 PPM	6		
	CBD / Cannabidiol	C21H30O2	12.739	12.640		313.2173	-0.57 PPM	2		
	CBN / Cannabinol	C21H26O2	12.282	12.173		309.1860	-0.08 PPM	2		
	Embelin	C17H26O4	6.280	6.191		293.1758	0.49 PPM	2		
i i	IsobutyIparaben	C11H14O3	3.574	3,497		193.0870	1.77 PPM	2		
	Thebaol	C16H14O3	6 165	6.095		253 0870	4 44 PPM	2		
	Diphenylmethoxyacetic acid	C15H14O3	5 558	5 458		241 0870	0.18 PPM	2		
	Fenoprofen	C15H14O3	5 558	5 453		241 0870	0.18 PPM	2		
	Nahumetone	C15H16O2	6 727	6 649		277 1078	-0.41 PPM	2		
	BPA / Bisphenol A	C15H16O2	6 727	6 6 2 9		227.1078	-0.41 PPM	2		
	Zingerone	C11H14O3	3 574	3 505		103 0870	1 77 DDM	2		
	Mosonin (Onianyl)	C10H1004	2 6 2 0	2 500		102 0506	0.20 DDM	2		
	Aspirin (Acobylealigylic acid)	C10/11004	4 520	1 127		170.0250	0.25 FFM	2		
	Coffoic acid	C9H6O4	4.520	4.437		179.0330	-0.20 PPM	2		
	Carreic aciu	C9H0U4	7.520	2,006		179.0330	-0.20 PPM	2		
	2 Dhanulahanal		3.962	3.900		1/7.06/0	0.79 PPM	2		
	2-Frienyphenor	C12H100	7.540	7.770		167.0009		2		
	Homogentisic acid		3.4/1	3.348		167.0350	-0.37 PPM	2		
	Metnyisalicyiate		2.691	2.595		151.0401	0.90 PPM	3		
+	Drimide	C18H24I3N3U8	2.179	0.393		789.8625	0.20 PPM	2		
	Primidone	C12H14N2O2	4.406	0.717		217.0983	4.33 PPM	1		
	5-Methylbenzotriazole	C/H/N3	3.853	0.536		132.0567	0.05 PPM		225 0240	
+	Estriol	C18H24O3	4.707	0.000		287.1653	-0.73 PPM	3	225.9218	
	Oxazepam	CI5HIICIN2O2	6.267	0.261		285.0436	-1.38 PPM	1		
	Isoproturon	C12H18N20	6.942	0.186		205.1346	-0.42 PPM	1		
	Diuron		0.517	0.250		231.0097	1.20 PPM	1		
	Napioxen Butul 4 hudrova honzonto (Butulnova honzo	C14H14O3	0.035	0.390		229.08/0	0.92 PPM	1		
	Ethinylestradial (EE2)	C11II1403	0.913	0.403		193.08/0	1.19 PPM	2	100 2002	
+	Ethinylestradiol (EE2)	C20H24O2	7.343	0.053		295.1/04	U.55 PPM	2	166.2993	
+	Estrone (E1)	C18H22U2	7.441	0.001		269.154/	-1.03 PPM	2	163.2462	
+	2-Phenyiphenol (Orthophenyiphenol)	C12H100	7.546	0.033		169.0659	0.00 PPM	2	215.1786	
+	Fluaioxonii	C12H6F2N2O2	7.937	0.238		247.0325	-0.16 PPM	3		
+	Gemtibrozil	C15H22O3	8.838	0.001		249.1496	0.15 PPM	3	382.1097	
	Merenamic acid	C15H15NO2	/.195	1./20		240.1030	0.15 PPM	3		
	Iriclosan	C12H7Cl3O2	10.151	0.536		286.9439	0.02 PPM	1		

図9.ターゲット化合物をスパイクした下水汚泥のネガティブイオン化モードで検出された推定上で同定された化合物

結論

このワークフローにより、下水サンプルにスパ イクしたターゲット化合物を定量できました。 分解能とダイナミックレンジが向上した 6546 LC/Q-TOF システムにより、PCDL からスク リーニングした 4,856 個の化合物から、8 個 の化合物がネガティブイオン化モードで高い 信頼性で確認され、同様に 18 個がポジティブ モードで確認されました。159 個のポジティブ イオン化化合物と 47 個のネガティブイオン化 化合物では、RT またはスペクトルを確認する 必要があります。確認が必要な化合物の多く が大麻の使用に関連していましたが、これは サンプリング場所での適法な使用です。キサン トフモールのようなポジティブおよびネガティ ブイオン化モードでスペクトルが精査されてい る一部の化合物が、ネガティブおよびポジティ ブイオン化モードの両方で検出されており、同 定の信頼性がさらに向上しています。チアベ ンダゾールなど、リテンションタイムが予測さ れた化合物も検出されました。リテンションタ イムの予測と新規汚染物質が掲載されたアジ レントの PCDL をさらに精査することで、存在 する可能性のある汚染物質の検出を引き続き 遡及的に実施できます。さらに多くの RT を予 測または見積もることにより、確認対象の化 合物の数が大幅に減少する可能性があります が、このワークフローでは SANTE によるサス ペクト化合物の同定基準を満たす可能性がな い多数の化合物を排除することにより、確認 の負担を減らしています。

環境毒性物質のリストは拡大しており、定量 能力への影響を最小限にして新しい化合物を 遡及的にモニタリングする機能を提供するこ とにより、ノンターゲット分析で従来の LC-TQ 分析手法全体の価値を高めています。

参考文献

- Black, G. P.; Anumol, T.; Young, T. M.; Analyzing a broader spectrum of endocrine active organic contaminants in sewage sludge with High Resolution LC-QTOF-MS suspect screening and QSAR toxicity prediction (In preparation).
- European Commission, Directorate General for Health and Food Safety. SANTE/11813/2017. 21–22 November 2017 rev.0
- Agilent SureMass Technical Overview, Agilent Technologies Technical Overview, publication number 5991-8048EN, 2017.
- Rennie, E. E.; *et al.* Agilent Technologies, publication number 5991-8580EN, **2017**.
- Berset, J. D.; Rennie, E. E.; Glauner, T. Agilent Technologies, publication number 5991-6627EN, 2016.

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. 2019 Printed in Japan, March 26, 2019 5994-0750JAJP

