

ICP-MS および統計モデリングを用いた 元素プロファイリングによる米の真正性判定

Agilent 7900 ICP-MS および Agilent Mass Profiler Professional ソフトウェアを用いた米の原産地の特定



はじめに

中国は世界最大の米生産国であり、米は世界人口のほぼ半数にとって重要な主食です。世界における 米の需要は非常に大きいため、高価値の品種は食品詐欺の標的となっており、金銭的利益のために高 価な食品への偽和物混入や偽ラベルの貼り付けが日常的に行われています。中国の米価は産地に大き く依存するため、生産者は地理的表示(GI)を利用して市場で製品を差別化します。原産地をセールス ポイントとする米で、生産可能な量を上回る需要がある場合は必ず、偽和物混入の被害に遭いやすくな ります。食品偽装の方法には、本物の製品を低品位の製品として部分的または完全に置き換える方法 や、不正確な製品ラベル付けを行う方法などがあります。食品偽装によって金銭的な影響を受ける可 能性があることから、農家、小売業者、消費者にとって、米の原産地の特定は重要です(1)。食品の産 地と真正性の検証は、食品安全性とコンプライアンスにおいても重要です。食品供給業者、製造業者、 小売業者が、汚染された食品や不正なラベルの付いた製品を供給していることが判明した場合、罰金 を科される可能性や起訴される可能性があります。

著者

Fei Xu¹, Fanzhou Kong¹, Hong Peng¹, Guangtao Zhang¹, and Shuofei Dong²

 Mars Global Food Safety Center, Beijing, China
 Agilent Technologies Co. Ltd (China), Beijing, China 食用作物と加工食品の元素組成は、植物の種類、土壌の種類、環境要因のほか、灌漑や肥料の使用といった農業慣行によって異なります。多くの場合、食品中の多岐にわたる元素の分析は、マイクロ波によるサンプルの酸分解の後に ICP-MS を用いて行われます (2-4)。その結果から各サンプルの「元素フィンガープリント」が得られ、それによってサンプルの原産地が示されます。この情報は真正性判定に有用で、茶や香辛料などの食品に適用されています (5、6)。

Agilent ICP-MS システムには、オクタポールリアクションシステム (ORS⁴) とコリジョン/リアクションセル (CRC) が搭載されています。こ れはヘリウム (He) コリジョンモードを使用して一般的な多原子イオン 干渉を完全に除去できるように最適化されています。ヘリウムモードで は、同じセル設定を多くの元素に使用できるため、食品の真正性判定の 研究に必要な高品質のデータセットが得られます。Agilent ICP-MS の データは、大量の MS データを処理するように設計された多変量解析ソ フトウェアパッケージである Agilent Mass Profiler Professional (MPP) と互換性があります。MPP にはデータ分類方法の選択肢が含まれてい るため、分析者は堅牢なモデルを構築して、未知のサンプルの属性(サ ンプルの産地など)を予測できます。

本研究では、原産地が判明している 90 種類の米サンプルの微量元素 データを、Agilent 7900 ICP-MS を使用して取得しまし、次に、大量の サンプルデータセットを MPP ソフトウェアによって処理しました。主成 分分析 (PCA)を使用して各サンプルの産地を区別可能と判断した後、 MPP ソフトウェアを用いて、さまざまなクラス予測アルゴリズムにより予 測モデルを構築しました。これらの予測モデルを使用して、24 個の「未 知」のテストサンプルの産地を特定しました。

実験方法

装置

すべての測定に、Agilent 7900 ICP-MS と Agilent SPS 4 オートサンプ ラを使用しました。7900 には、ガラス製同軸ネブライザ、石英製ダブル パススプレーチャンバ、内径 2.5 mm インジェクタ付き石英トーチ、Ni イ ンタフェースコーン、および ORS⁴ セルが含まれています。装置の制御に は Agilent ICP-MS MassHunter ソフトウェアを使用し、データ解析は MPP ソフトウェアを用いて実行しました。ここに記載した機能は、MPP バージョン 15.0 以降で使用できます。主な装置操作パラメータを表 1 に 示します。 表1. Agilent 7900 ICP-MS の使用条件

パラメータ	設定値
RF 出力(W)	1550
サンプリング深さ(mm)	8
ネブライザガス (L/min)	1.16
レンズチューン	オートチューン
He 流量(mL/min)	5
KED (V)	5

検量用標準溶液

24 元素の検量用標準溶液は、希釈に5% (v/v) 硝酸 (HNO₃) 溶液 を使用して Agilent 標準溶液から調製しました。含まれる標準は、混 合標準溶液-2A (部品番号 8500-6940)、混合標準溶液-4 (部品番号 8500-6942)、および環境検量用標準溶液(部品番号 5183-4688) で す。ロジウム (Rh) 内部標準 (ISTD) 溶液は、単一元素の Rh 標準(部 品番号 5190-8509) から調製しました。水銀 (Hg) を除く検量用標準 溶液は、分析前の 24 時間以内に事前混合しました。Hg は使用直前に 混合溶液に添加しました。

試薬、参照物質、サンプル

硝酸 (69 %) は、Merck Millipore (ドイツダルムシュタット) から購入しました。脱イオン水 (DIW、18.2M Ω cm) は、Milli-Q システム (Millipore、米国マサチューセッツ州) から入手しました。

米粉標準物質 (SRM) 1568b は、米国国立標準技術研究所 (NIST、 米国メリーランド州ゲーサーズバーグ)から購入しました。産地が判明し ている 90 の米サンプル (GI 指定製品)は、中国の5つの省の米加工工 場から直接収集しました。遼寧省から収集した2種類の30サンプルを 除いて、各省から15サンプルを集めました。収集した場所は、黒竜江省 五常市、遼寧省盤錦市 (異なるタイプのサンプル2ロット)、江蘇省射陽 県、湖北省京山市、広西チワン族自治区貴港市でした。サンプルには、そ れぞれ HLJ、LN-1、LN-2、JS、HB、GX のラベルを付けました。

サンプル前処理

分解用容器を含むすべてのラボ材料を30% (v/v) HNO₃ 溶液に24時 間浸し、使用前に DIW で3回すすぎました。米サンプル約0.5gを計量 してポリテトラフルオロエチレン (PTFE) 分解用容器に移し、6 mL の HNO₃ と混合しました。容器をドラフト内に一晩置いて、サンプルを事前 分解しました。次に、酸分解のためにマイクロ波オーブン (Anton Paar、 オーストリア) に移しました。酸分解には、表2に示す加熱プログラムを 使用しました。 分解後、溶液を室温まで冷却し、DIW で 50 mL に希釈しました。分解メ ソッドを検証するために NIST-1568b SRM を米サンプルのすべてのバッ チで分析しました。

表 2. マイクロ波分解の代表的なパラメータ

ステージ	時間	温度
昇温	15 分	室温から 180 ℃
保持	20 分	180 °C
冷却	約 45 分	180 ℃から室温

結果と考察

検出限界 (DL) およびバックグラウンド相当濃度 (BEC)

24 元素の ICP-MS MassHunter キャリブレーションから計算した 7900 ICP-MS の代表的な DL および BEC を表 3 に示します。

表 3. 本研究で測定した 24 元素の DL と BEC

元素	DL (µg/L)	BEC (µg/L)	元素	DL (µg/L)	BEC (µg/L)
11 B	0.33	0.064	60 Ni	0.012	0.047
23 Na	0.091	1.03	63 Cu	0.0011	0.016
24 Mg	0.092	0.075	66 Zn	0.018	0.18
27 Al	0.37	0.47	75 As	0.0020	0.0020
39 K	0.39	11.90	78 Se	0.048	0.019
44 Ca	0.030	0.137	85 Rb	0.00001	0.0009
47 Ti	0.0064	0.0021	88 Sr	0.0015	0.0006
51 V	0.0015	0.0007	95 Mo	0.0011	0.0019
52 Cr	0.0021	0.014	111 Cd	0.0039	0.0049
55 Mn	0.0037	0.0042	137 Ba	0.0033	0.0069
56 Fe	0.022	0.13	201 Hg	0.0001	0.0009
59 Co	0.0005	0.0062	208 Pb	0.0097	0.12

SRM の回収率

NIST-1568b 米粉 SRM を各バッチランで 2 ~ 3 回分析し、メソッドの正 確度と精度を検証しました。4 回のバッチランの結果を表 4 に示します。 平均濃度は、Hg を除くすべての元素の認定濃度(回収率 80 ~ 120 %) とよく一致しました。Hg の回収率が低いのは、5 ~ 8 % HNO₃ 中の Hg が低濃度で不安定であったためと考えられます(7)。

サンプルの分解と安定化に使用される酸混合物に HCI を添加すると、 Hg の回収率が大幅に向上する可能性があります。HCI により、化学的 に不安定な Hg などの元素を溶液中に確実に保持できます。形成された CI ベースの多原子イオン干渉は、Agilent ICP-MS システムで標準の He セルモードを使用して簡単に制御できます(8)。

- +		مر بر الله من المراجع ا	1	
元素	認定濃度 (µg/g)	平均測定濃度	標準偏差	平均回収率
		(µg/g)	(µg/g)	(%)
24 Mg	559 ± 10	540	27	97
27 Al	4.21 ± 0.34	4.11	0.15	98
39 K	1282 ± 11	1276	62	100
44 Ca	118.4 ± 3.1	114.3	8.3	97
55 Mn	19.2 ± 1.8	18.3	1.0	95
56 Fe	7.42 ± 0.44	7.30	0.30	98
59 Co*	0.0177 ± 0.0005	0.0170	0.0010	96
63 Cu	2.35 ± 0.16	2.29	0.15	98
66 Zn	19.42 ± 0.26	17.73	1.81	91
75 As	0.285 ± 0.014	0.269	0.020	94
78 Se	0.365 ± 0.029	0.351	0.026	96
85 Rb	6.198 ± 0.026	5.352	0.211	86
95 Mo	1.451 ± 0.048	1.405	0.110	97
111 Cd	0.0224 ± 0.0013	0.0197	0.0022	88
201 Hg	0.00591 ± 0.00036	0.00395	0.00127	67
208 Pb*	0.008 ± 0.003	0.009	0.001	116

表 4. NIST-1568b 米粉 SRM 中の認定元素の回収率と再現性データ、n = 10

*参考値:完全に検証された値ではありませんが、認証値と互換的に使用できます。

ISTD 安定性

Rh は、ISTD として使用しました。これは、Sc や Ga などの他の ISTD 元素と異なり、Rh は米サンプル中に存在するとは予想されなかったた めです。Rh 標準溶液(1 mg/L)は、T 型コネクタを使用してサンプルと 混合しました。すべての米サンプルの半分からなる分析シーケンスを、10 時間にわたって繰り返し分析しました。図 1 に赤い点線で示すように、す べての ISTD の回収率は ± 20 % の範囲内でした。この結果は、食品分 解物の長期分析を必要とする真正性研究において、7900 ICP-MS が堅 牢性と高いマトリックス耐性を備えていることを示しています。



図 1. Agilent 7900 ICP-MS を使用した 10 時間にわたるサンプル分析の ISTD 回収率

MPP ソフトウェアを用いたデータマイニング

多元素データバッチファイル (米サンプル 90 個、元素 24 種類) を組み 合わせて、統計解析のために MPP ソフトウェアにインポートしました。こ のソフトウェアには、t 検定などのさまざまな検定、分散解析 (ANOVA)、 モデル構築アルゴリズム、相関解析、クラスタリング解析ほか、R アルゴ リズムの拡張機能が含まれています。PCA (教師なし手法)を実行して、 サンプルグループ間に有意差があるかどうかを検証しました。PCA 手法 は、グループの分離に対する化合物の相対的な寄与を評価します。3 次元 (3D) PCA スコアを図 2 に示します。

分散比の合計の65%を、最初の3つの次元、PC1、PC2、およびPC3 が占めました(図2)。24元素を使用して、5つの産地の90の米サンプ ルが適切な程度で分離されました。HB、GX、およびLN-1のサンプルは、 他のサンプルから明確に分離されましたが、LN-2とHLJのサンプルに は多少の重なりがありました。MPP ソフトウェアによる PCA ローディン グプロットに基づくと、違いは主にPC1のB、Na、Zn、Cdと、PC2の Al、Fe、Mn、Sr、およびPC3のNa、Cu、Se、Rb に起因するものでした。 この結果は、元素濃度データを使用した産地予測モデルの構築の実現可 能性を示しています。これにより、未知の米サンプルの原産地の予測が可 能になると考えられます。



図 2.5 つの場所 (遼寧省の2つの異なる品種を含む)からの 90 の米サンプルの 3D-PCA スコア。グラフの軸は、PCA 結果の上位3つの成分を表します。

クラス予測解析

クラス予測解析は、未知の新しいサンプルを以前に決定されたグループ に偏りなく割り当てることができる有用な手法です。ワインやビールなど の複雑なサンプルの品質管理において、ますます価値のあるツールにな りつつあります (9)。

MPP には、複数のクラス予測アルゴリズムが含まれています(図3)。 アルゴリズムの選択肢には、SIMCA (Soft Independent Modeling of Class Analogy)、ランダムフォレスト、線形判別分析(LDA)、PLS 判別 分析(PLSDA)、決定木、サポートベクターマシン(SVM)、単純ベイズ、 およびニューラルネットワークなどがあります。分類モデルの構築には、 これらのクラス予測アルゴリズムのいずれかを使用できます。



図3. クラス予測アルゴリズムの選択肢が表示された Agilent MPP ソフトウェア画面

MPP での予測モデルの構築

5 つの各場所 (遼寧省からは 2 セット) からの 15 の米サンプルの うち11 個をランダムに選択し、4 つの異なるクラス予測アルゴリズム (PLSDA、SVM、LDA、SIMCA)を用いて予測モデルを構築しました。 各場所からの残り 4 つの米サンプル (合計 24 サンプル) は、予測モデル の精度をテストするための「未知」として使用しました。MPP では、6 ス テップの自動ワークフローに従うだけで予測モデルを構築でき、不要な ステップはソフトウェアが自動的にスキップします。

予測モデルのテスト

PLSDA、SVM、および LDA モデルを使用して得られた予測産地の結果 を表 5 に示します。24 サンプルすべての産地が、3 つの予測モデルすべ てで正しく特定されました。表 5 の値が大きい(陰影が濃い)ほど、正解 との一致の信頼度が高いことを示しています。

		予測モデル							
サンプル ID 実際の産地		PLSDA		SVM		LDA			
		予測された産地	信頼度	予測された産地	信頼度	予測された産地	信頼度		
GX-101	広西チワン族自治区貴港市	広西チワン族自治区貴港市	0.762	広西チワン族自治区貴港市	0.742	広西チワン族自治区貴港市	0.776		
GX-102	広西チワン族自治区貴港市	広西チワン族自治区貴港市	0.476	広西チワン族自治区貴港市	0.758	広西チワン族自治区貴港市	0.658		
GX-103	広西チワン族自治区貴港市	広西チワン族自治区貴港市	0.606	広西チワン族自治区貴港市	0.804	広西チワン族自治区貴港市	0.830		
GX-104	広西チワン族自治区貴港市	広西チワン族自治区貴港市	0.769	広西チワン族自治区貴港市	0.769	広西チワン族自治区貴港市	0.865		
HB-101	湖北省京山市	湖北省京山市	0.632	湖北省京山市	0.774	湖北省京山市	0.667		
HB-102	湖北省京山市	湖北省京山市	0.739	湖北省京山市	0.797	湖北省京山市	0.760		
HB-103	湖北省京山市	湖北省京山市	0.770	湖北省京山市	0.727	湖北省京山市	0.811		
HB-104	湖北省京山市	湖北省京山市	0.531	湖北省京山市	0.757	湖北省京山市	0.850		
HLJ-101	黒竜江省五常市	黒竜江省五常市	0.423	黒竜江省五常市	0.415	黒竜江省五常市	0.733		
HLJ-102	黒竜江省五常市	黒竜江省五常市	0.651	黒竜江省五常市	0.731	黒竜江省五常市	0.901		
HLJ-103	黒竜江省五常市	黒竜江省五常市	0.659	黒竜江省五常市	0.777	黒竜江省五常市	0.873		
HLJ-104	黒竜江省五常市	黒竜江省五常市	0.635	黒竜江省五常市	0.724	黒竜江省五常市	0.802		
JS-101	江蘇省射陽県	江蘇省射陽県	0.637	江蘇省射陽県	0.766	江蘇省射陽県	0.834		
JS-102	江蘇省射陽県	江蘇省射陽県	0.494	江蘇省射陽県	0.625	江蘇省射陽県	0.866		
JS-103	江蘇省射陽県	江蘇省射陽県	0.526	江蘇省射陽県	0.735	江蘇省射陽県	0.918		
JS-104	江蘇省射陽県	江蘇省射陽県	0.562	江蘇省射陽県	0.824	江蘇省射陽県	0.914		
LN1-101	遼寧省盤錦市 -1	遼寧省盤錦市 -1	0.576	遼寧省盤錦市 -1	0.711	遼寧省盤錦市 -1	0.881		
LN1-102	遼寧省盤錦市 -1	遼寧省盤錦市 -1	0.738	遼寧省盤錦市 -1	0.743	遼寧省盤錦市 -1	0.893		
LN1-103	遼寧省盤錦市 -1	遼寧省盤錦市 -1	0.546	遼寧省盤錦市 -1	0.744	遼寧省盤錦市 -1	0.450		
LN1-104	遼寧省盤錦市-1	遼寧省盤錦市-1	0.594	遼寧省盤錦市 -1	0.816	遼寧省盤錦市 -1	0.838		
LN2-101	遼寧省盤錦市 -2	遼寧省盤錦市-2	0.769	遼寧省盤錦市 -2	0.717	遼寧省盤錦市 -2	0.832		
LN2-102	遼寧省盤錦市 -2	遼寧省盤錦市-2	0.593	遼寧省盤錦市 -2	0.731	遼寧省盤錦市-2	0.774		
LN2-103	遼寧省盤錦市 -2	遼寧省盤錦市-2	0.763	遼寧省盤錦市 -2	0.734	遼寧省盤錦市 -2	0.863		
LN2-104	遼寧省盤錦市 -2	遼寧省盤錦市-2	0.535	遼寧省盤錦市 -2	0.727	遼寧省盤錦市 -2	0.745		

表 5.24の「未知」米サンプルの産地予測結果のまとめ

SIMCA モデルを使用して得られた、同じテストサンプルの結果を表 6 に 示します。SIMCA モデルは、「距離」を使用して結果を報告します。値が 小さいほど、サンプルが分類されたグループに属する可能性が高くなりま す。SIMCA モデルの予測は、1 つの米サンプル(HLJ-101)を除いて、 すべて正しいものでした。HLJ-101 は LN-2 グループの可能性が最も高 いという予測となり、PLSDA および SVM モデルでも低い信頼値となり ました(表 5)。PCA の結果は、HLJ と LN-2 の元素プロファイルにいく つかの類似点があることを示唆しています(図 2)。

サンプル ID	実際の産地	予測アルゴリズム					
		SIMCA					
		距離(広西チワン族 自治区貴港市)	距離 (黒竜江省五常市)	距離 (湖北省京山市)	距離 (江蘇省射陽県)	距離 (遼寧省盤錦市 -1)	距離 (遼寧省盤錦市 -2)
GX-101	広西チワン族自治区貴港市	0.857	8.050	2.901	8.363	7.635	3.254
GX-102	広西チワン族自治区貴港市	0.203	9.422	4.627	7.672	8.797	4.865
GX-103	広西チワン族自治区貴港市	0.221	11.803	2.816	8.288	8.953	3.522
GX-104	広西チワン族自治区貴港市	0.195	8.879	3.965	7.573	8.219	3.801
HB-101	湖北省京山市	3.759	11.049	0.255	9.807	6.379	4.664
HB-102	湖北省京山市	5.279	10.986	0.527	10.328	5.746	5.411
HB-103	湖北省京山市	5.179	10.519	0.343	9.809	5.585	5.048
HB-104	湖北省京山市	4.126	10.093	0.367	9.026	5.288	4.650
HLJ-101	黒竜江省五常市	4.806	3.262	4.684	4.202	3.043	1.841
HLJ-102	黒竜江省五常市	7.328	0.370	6.634	2.685	3.246	1.511
HLJ-103	黒竜江省五常市	8.362	0.407	7.177	2.515	3.702	1.475
HLJ-104	黒竜江省五常市	8.740	0.630	7.133	2.351	3.012	1.895
JS-101	江蘇省射陽県	14.174	6.839	16.351	0.141	2.901	4.514
JS-102	江蘇省射陽県	12.018	6.895	14.477	0.647	2.830	4.234
JS-103	江蘇省射陽県	12.831	6.750	16.327	0.292	3.003	2.786
JS-104	江蘇省射陽県	12.003	7.809	17.050	0.146	4.089	3.750
LN1-101	遼寧省盤錦市 -1	15.744	10.073	10.605	3.314	0.141	5.148
LN1-102	遼寧省盤錦市 -1	18.586	8.118	12.548	3.465	0.302	6.626
LN1-103	遼寧省盤錦市 -1	24.607	12.810	22.225	3.157	0.892	7.842
LN1-104	遼寧省盤錦市 -1	17.242	12.822	12.874	5.091	0.193	6.848
LN2-101	遼寧省盤錦市 -2	11.641	2.939	8.208	2.757	2.114	0.400
LN2-102	遼寧省盤錦市 -2	11.823	2.944	9.358	3.014	2.573	0.266
LN2-103	遼寧省盤錦市 -2	8.010	2.334	8.650	2.498	3.692	0.379
LN2-104	遼寧省盤錦市 -2	9.582	3.681	9.425	2.805	3.444	0.570

表 6. SIMCA モデルによる 24 の「未知」の米サンプルの産地予測の結果。緑:正しい予測、茶色:誤った予測、赤:実際の産地

結論

ヘリウムコリジョンモードで動作する ORS⁴ セルを備えた Agilent 7900 ICP-MS により、食品の真正性分析に不可欠な品質、堅牢性、安定性に 優れたデータが得られることを示しました。すべての分析対象物にシング ルセルガスモードを使用することで、90 個の米サンプル分解物中の 24 種類の元素を検出することができました。90 個の米サンプルのうち 66 個のデータを使用して、残り 24 サンプルの原産地を特徴付ける予測モデ ルを構築しました。

大量のデータセットの処理には、Agilent MPP ソフトウェアを使用しました。PCA を使用したデータ解析では、米の元素組成が原産地の影響を 受けており、5つの地理的領域をかなり的確に区別できることが示されま した。

MPP ソフトウェアの簡単な手順に従って、4 つの予測モデルを作成しま した。モデルのトレーニングには ICP-MS のデータを使用し、「未知」の サンプルを使用してテストしました。本研究では、単一の予測モデルでは なく複数の予測モデルを使用することで、24 個の「未知」のサンプルす べての産地を正しく予測できることが示されました。複数モデルによるア プローチは、より正確な予測結果とサンプルの元素プロファイルのより良 い理解につながる可能性があります。

本研究で紹介した元素プロファイリングメソッドは、米やその他の高価値 食品の原産地を特徴づけて、食品真正性のルーチン分析を実現する可能 性を有しています。

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. 2021 Printed in Japan, September 6, 2021 5994-4043JAJP DE44425.119375

参考文献

- 1. Dion M.A.M.Luykx, Saskia M. van Ruth, An overview of analytical methods for determining the geographical origin of food products, *Food Chemistry*, **2008**, 107, 897–911
- 2. Jenny Nelson, Elaine Hasty, Leanne Anderson, Macy Harris, US FDA EAM 4.7 ICP-MS メソッドに従った食品中の元素測定, Agilent publication, <u>5994-2839JAJP</u>
- Shuofei Dong, Jenny Nelson, Michiko Yamanaka, Routine Analysis of Fortified Foods using the Agilent 7800 ICP-MS, Agilent publication, <u>5994-0842EN</u>
- Kazuhiro Sakai, Junichi Takahashi and Ed McCurdy, 自動メソッド作成機能を備えた Agilent 7900 ICP-MS の食品中 ルーチン微量金属元素分析への応用, Agilent publication, <u>5991-</u> <u>4556JAJP</u>
- Jenny Nelson, Helene Hopfer, Authentication of Specialty Teas: An Application Note.*Food Qual.Safety*, **2018**, December, accessed February 2021, <u>https://www.foodqualityandsafety.</u> <u>com/article/specialty-tea-authentication/</u>
- Jenny Nelson et al., Determining the metal content of the spices and identifying the country of origin, *Food Qual. Safety*, **2019**, May, accessed February 2021, <u>https://www. foodqualityandsafety.com/article/authentic-spicesidentifying-country-origin/
 </u>
- U.S. EPA.2014.Method 6020B (SW-846): Inductively Coupled Plasma-Mass Spectrometry, Revision 2.Washington, DC.<u>https://www.epa.gov/sites/production/files/2015-12/</u> <u>documents/6020b.pdf</u>
- 8. 効率に優れた水銀の低濃度分析, Agilent publication, <u>5990-</u> <u>7173JAJP</u>
- Spiros A. Drivelos, Constantinos A Georgiou, Multielement and multi-isotope-ratio analysis to determine the geographical origin of foods in the European Union, *Trends Anal.Chem*, **2012**, 40, 38–51

