

マイクロ波プラズマ原子発光分光 によるエタノール燃料中の Cr、Ni、Pb、V の測定

アプリケーションノート

エネルギー・燃料

著者

George L. Donati*, Renata S. Amais*, Daniela Schiavo[†], Joaquim A. Nóbrega*

* Group of Applied Instrumental Analysis, Departament of Chemistry, Federal University of São Carlos, São Carlos, SP, Brazil

† Agilent Technologies São Paulo, SP, Brazil



概要

マイクロ波プラズマ原子発光分光分析装置 (MP-AES) による簡単な希釈手順を用いた、エタノール燃料中 Cr、Ni、Pb、Vの測定方法を紹介します。サンプル前処理が簡単で、特殊なガスや高価なガスは必要ありません。この分析で得られる検出下限 (0.3~40 µg/L) はいずれも、燃料燃焼による環境負荷やエンジン性能における燃料汚染の影響に関する主要法規制の要件を満たしています。



はじめに

エタノール燃料は、1970年代の最初の石油危機以来、自動車で 広く使用されるようになっています。最近では、数多くのフレッ クス燃料エンジンの発売と、再生可能エネルギーの使用による 環境上の利点により、世界中でエタノールの生産と消費が拡大 しています。ブラジルでは、エタノール燃料はサトウキビからつ くられます。サトウキビは収穫量が多く、持続可能なエネルギー の1つに数えられます[1]。

燃料中に金属が存在すると、酸化分解反応により、エンジン 性能が低下したり、燃料品質が悪化したりすることがありま す[2]。また、サトウキビの生育した土壌の分解により、一部の 毒性元素が自然にエタノール中に混入することもあります。あ るいは、製造、保管、輸送の際に毒性元素が燃料に入りこむケー スもあります。そのため、こうした毒性元素により、燃料燃焼後 の空気汚染が拡大するおそれがあります[3]。

このアプリケーションノートでは、Agilent 4100 マイクロ波プラ ズマ原子発光分光分析装置 (MP-AES) を用いたエタノール燃料 の直接分析による Cr、Ni、Pb、V の測定方法を紹介します。この 装置は、磁気的に結合したマイクロ波エネルギーにより生じる 窒素プラズマをベースにしています。そのおもな利点としては、 ランニングコストやメンテナンスコストの削減が挙げられます。 窒素ガスジェネレータとエアコンプレッサだけで十分に機器を 動作できるため、他のガスは必要ありません。この実験では、 1 % v/v HNO₃ 水溶液を用いたエタノールの希釈により、正確な 分析結果を簡単に得ることができました。

実験手法

使用機器

Agilent 4100 MP-AES を使用しました。サンプル導入システムの 構成は、溶媒耐性チューブ、ダブルパスサイクロニックスプレー チャンバ、不活性 OneNeb ネブライザです。水溶液による希釈 以外の前処理をおこなわずに、燃料サンプルを直接導入するた め、外部ガスコントロールモジュール (EGCM)を用いて空気を 窒素プラズマに導入し、トーチおよびプレ光学ウィンドウでの 炭素析出を防止しました。空気を導入することで、プラズマを 安定化させ、バックグラウンド発光を低減する効果も得られ ます。 Agilent MP Expert ソフトウェアのオートバックグラウンド補正 機能を用いて、バックグラウンド補正を自動的におこないまし た。この機能は、各元素のバックグラウンドスペクトルを記録お よび保存し、差し引きます。バックグラウンドスペクトルはブラ ンク溶液から得られるものです。その後、すべてのサンプル溶 液から、このスペクトルが差し引かれます。また、モニタリング する各波長について、ネブライザ圧力や観測ポジションなどの パラメータを自動的に最適化することもできます。表1と2に、 エタノール燃料サンプル中の Cr、Ni、Pb、V の測定に用いた機 器動作条件と設定を示しています。

表 1. エタノール燃料サンプルの直接分析に用いた Agilent 4100 MP-AES 動作条件

機器パラメータ	動作条件
ネブライザ	不活性 OneNeb
スプレーチャンバー	サイクロニックダブルパス
読み取り時間 (秒)	5
反復回数	3
安定化時間 (秒)	15
バックグラウンド補正	Auto

表 2. Cr、Ni、Pb、V の測定に用いたネブライザ圧力と EGCM 設定

元素	波長 (nm)	ネブライザ圧力 (kPa)	EGCM
Cr	425.433	240	高
Ni	352.454	180	高
Pb	405.781	100	高
V	437.923	240	高

試薬と標準溶液

すべての溶液の前処理には、サブボイリング蒸留システム (Milestone、ソリーゾレ、イタリア) によりあらかじめ精製 した硝酸 (Merck、ダルムシュタット、ドイツ) を使用しまし た。それぞれ 1000 mg/L の Cr、Ni、Pb、V を含む単元素原液 (Tec-Lab、Hexis、サンパウロ、SP、ブラジル)を用いて標準溶液を 作成し、添加実験をおこないました。分析グレードのエタノール (J. T. Baker、ピッツバーグ、NJ、USA)を用いて、検量線の 作成に使用する標準溶液のマトリックスマッチングをおこない ました。

サンプルとサンプル前処理

エタノール燃料サンプル (水和エタノール) をサン・カルロス (SP、ブラジル)のガソリンスタンドで入手しました。ブラジルの 法規則の規定では、水和エタノール燃料に含まれる水の最大許 容量は 4.9 % v/v です [4]。サンプルを HNO31% v/v で 10 倍に 希釈しました。1% v/v HNO3中にそれぞれ Cr、Ni、Pb、V を含む 有機標準溶液の適量を希釈し、外部キャリブレーションメソッ ドに使用する標準溶液を作成しました。各標準溶液にエタノー ルを加え、最終濃度を 10% v/v としました。

結果と考察

性能指標

サンプル分析前に、各分析対象物の分析性能を評価しました。 バックグラウンド相当濃度 (BEC)、シグナル/バックグラウンド 比 (SBR)、各 10 回の連続ブランク測定の相対標準偏差から、 検出下限 (LOD) と定量下限 (LOQ) を算出しました。表 3 に、得 られた LOD および LOQ を示しています。これらのデータか らは、フレーム原子吸光分光法 (FAAS) などと比べた場合の 4100 MP-AES の優れた分析性能が見てとれます。マイクロ波プ ラズマは、特に Cr や V などの難揮発性元素で大きな効果を発 揮します。このケースでは、プラズマでの到達温度が高いため、 優れた LOD が得られます。亜酸化窒素-アセチレンフレームを 用いる FAAS 分析のように特殊なガスを導入する必要はありま せん [5]。

エタノール燃料サンプルの直接分析の精度を評価するために、 添加実験をおこないました。結果を表4に示しています。分析 したすべての元素で、回収率は92~108%でした。このことは、 有機化合物やCu、Na、Feなどの元素により生じるマトリック ス効果が小さいことを示しています。 **表 3**. MP-AES によるエタノール燃料中 Cr、Ni、Pb、V 測定の性能指標 * 機器の検出下限と定量下限

+ サンプル希釈 (HNO₃1% v/v 中に 1:9 v/v エタノール燃料) を考慮し た検出下限

元素	LOD* (µg/L)	LOQ* (µg/L)	サンプルの LOD [†] (µg/kg)
Cr	0.7	2.2	9
Ni	16	52	200
Pb	40	130	490
v	0.3	0.9	4

表4.エタノール燃料サンプル中Cr、Ni、Pb、V測定の添加実験

元素	添加 (µg/L)	検出 (μg/L)	回収率 (%)
Cr	20	21.2 ± 1.2	106
	100	95.1 ± 1.2	95
	500	460 ± 30	92
Ni	100	95.3 ± 0.8	95
Pb	400	430 ± 10	108
	1000	990 ± 10	99
V	20	19.8 ± 1.6	99
	100	98.4 ± 1.4	98
	500	460 ± 20	92

結論

Agilent 4100 MP-AES を用いたエタノール燃料の直接分析は、 ルーチン分析に容易に導入できるシンプルで効果的なメソッド です。3 種類のサンプルを分析したところ、いずれも分析対象 元素による汚染は見られませんでした(すなわち、元素濃度が 検出下限未満)。1 % v/v HNO3 を用いた簡単な希釈注入手順に より、Cr、Ni、Pb、V を正確に測定できました。EGCM により、トー チとプレ光学コンポーネントでの炭素析出が防止され、バック グラウンドシグナルの低減および精度向上という効果が得られ ました。

参考文献

- Silva, J. E., Silva, F. A., Pimentel, M. F., Honorato, R. S., Silva, V. L., Montenegro, M. C. B. S. M. & Araújo, A. N. (2006). A flow-batch internal standard procedure for iron determination in hydrated ethanol fuel by flame atomic absorption *spectrometry. Talanta*, *70*, 522–526.
- [2] Oliveira, M. F., Saczk, A. A., Okumura, L. L., Fernandes, A. P., Moraes, M. & Stradiotto, N. R. (2004). Simultaneous determination of zinc, copper, lead, and cadmium in fuel ethanol by anodic stripping voltammetry using a glassy carbon-mercury film electrode. *Anal. Bioanal. Chem., 380*, 135–140.
- [3] Saint'Pierre, T. D., Maranhão, T. A., Frescura, V. L. A. & Curtius, A. J. (2005). The development of a method for the determination of trace elements in fuel alcohol by electrothermal vaporization - inductively coupled plasma mass spectrometry using external calibration. *Spectrochim. Acta Part B, 60*, 605–613.
- [4] Brazilian National Agency of Petroleum, Natural Gas and Biofuels. Resolution ANP 7, D.O.U. 10.2.2011.
- [5] Amorim, F. A. C., Welz, B., Costa, A. C. S., Lepri, F. G., Vale, M. G. R. & Ferreira, S. L. C. (2007). Determination of vanadium in petroleum and petroleum products using atomic spectrometric techniques. *Talanta*, *72*, 349–359.

www.agilent.com/chem/jp

アジレントは、本文書に誤りが発見された場合、また、本文書の 使用により付随的または間接的に生じる損害について一切免責 とさせていただきます。

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。著作権法で許されている場合を除き、書面に よる事前の許可なく、本文書を複製、翻案、翻訳することは禁じ られています。

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2012 Published August 9, 2012 Publication number: 5991-0771JAJP

