

LC/MS によるリチウムイオン電池電解液の 品質管理



著者

Florian Rieck Agilent Technologies, Inc.

概要

リチウムイオン電池の効率と寿命は、電解液の品質に大きく依存します。現在使用されているほとんど のリチウムイオン電池の電解液成分には、水との反応によって分解しやすい性質があります。このアプ リケーションノートでは、頻繁に使用される電解液溶媒とその分解生成物の一部を分離して定量する HPLC-MS メソッドを紹介します。Agilent 1260 Infinity II ハイブリッドマルチサンプラを Feed 注入モー ドで使用した場合、サンプルが強溶媒であるにもかかわらず、ピークの歪みは確認されませんでした。 主要成分の検出下限は 30 ppm 未満、分解生成物の検出下限は 5 ppb 未満でした。

はじめに

リチウムイオン電池 (LIB) は、一般消費者向けのポータブル電子機器や 電気自動車で広く使用されています。より効率的で軽量な大容量電池の 開発は、大きな研究分野です。LIB の効率に関係する重要な要素は、電 解液とその成分です。電解液は、有機炭酸塩溶媒に溶解したリチウム塩 (最も一般的なものは LiPF₆)で構成されています。¹ これらの成分は水 分の影響を受けやすく、高温で充電サイクル数が多くなると経時的に分解 します。したがって、LIB の長期的な効率を判断するためには、主要成分 と分解生成物の測定が重要になります。

エチレンカーボネート (EC) とジアルキルカーボネートの混合物が電 解液中に存在する場合、必ずオリゴマーカーボネートに反応します (図 1)。EC、ジメチルカーボネート (DMC)、ジエチルカーボネート (DEC) を含む電解液は、例えば、ジメチル 2,5-ジオキサへキサンジオエー ト (DMDOHC、R = CH₃) とジエチル 2,5-ジオキサへキサンジオエート (DEDOHC、R = C₂H₅) に分解されます。



図 1. EC およびジアルキルカーボネートからのオリゴマーカーボネートの生成。 参考文献 4 から改変しています。

このアプリケーションノートでは、一般的な電解液溶媒とその分解生成物 を分離して定量する HPLC-MS メソッドを紹介します。この分離メソッド には、EC、DMC、DEC、DMDOHC、DEDOHC の他に、同じく一般的な 電解液溶媒である、エチルメチルカーボネート (EMC) やプロピレンカー ボネート (PC) も含まれています。リチウム塩中の元素不純物の測定に ついては、参考文献 2 および 3 を参照してください。

実験方法

使用装置

1260 Infinity II LC は、次のモジュールで構成しました。

- Agilent 1260 Infinity II フレキシブルポンプ (G7104C)
- Agilent 1260 Infinity II ハイブリッドマルチサンプラ (G7167C)
- Agilent 1260 Infinity II マルチカラムサーモスタット (G7116A)
- Agilent 1260 Infinity II ダイオードアレイ検出器 WR (G7115A)
- Agilent InfinityLab LC/MSD iQ (G6160A)

カラム

Agilent InfinityLab Poroshell 120 EC-C18、2.1 × 250 mm、4 μm (部 品番号 650750-902T)

ソフトウェア

Agilent OpenLab CDS リビジョン 2.6 以降

溶媒

移動相とサンプル溶媒として、Agilent InfinityLab アセトニトリル (ACN) LC/MS 用 (部品番号 5191-5101*)を使用しました。超純水 は、0.22 µm メンブレンユースポイントカートリッジ (Millipak)を装着し た Milli-Q Integral システムで製造しました。LC/MS グレードのギ酸は、 VWR (ダルムシュタット、ドイツ) から購入しました。

サンプル

参照電解液成分のエチレンカーボネート(EC)>99%、プロピレンカー ボネート(PC)99.7%、ジメチルカーボネート(DMC)>99%、エチル メチルカーボネート(EMC)99.9%、ジメチル2,5-ジオキサへキサンジ オエート(DMDOHC)98%、ジエチルカーボネート(DEC)>99%、ジ エチル2,5-ジオキサへキサンジオエート(DEDOHC)98%、およびジプ ロピルカーボネート(DPC)99%は、Merck(ダルムシュタット、ドイツ) から購入しました。EC、DMC、DEC(1:1:1、v:v)中の電池グレードの 1 M ヘキサフルオロリン酸リチウム溶液もMerckから入手し、電解液サ ンプルとして使用しました。

EC、PC、DMDOHC、DEDOHC については 10 ~ 10,000 ppb の範囲で、 DEC、DMC、EMC については 10 ~ 100,000 ppm の範囲で、すべての 参照電解液の検量線を作成しました。内部標準には DPC を使用しました。

技術的な理由から、サンプルと標準を、不活性ガス下のグローブボックス 内で取り扱うことはできませんでした。ただし、この制限は分解生成物の 生成に有利であることが判明しました。その検出がこのアプリケーション ノートの目的でした。

^{*}特定の国でのみ入手可能です。

メソッドの設定

表1. クロマトグラフィー条件

パラメータ	設定値				
移動相	A) 0.1 % ギ酸(FA)水溶液 B) 0.1 % ギ酸アセトニトリル溶液				
流量	0.4 mL/min				
グラジエント	0分 2% B 2分 2% B 11分 60% B 12分 95% B 14分 95% B 14.5分 2% B				
ストップタイム	20分				
注入量	1 μL				
Feed 注入	フィード速度:20 μL/min オーバーフィード容量:3.5 μL フィード溶媒 (S2):水+2% ACN+0.1% FA 内側洗浄:150 μL S1 再コンディショニング:S2				
ニードル洗浄(S1)	水:アセトニトリル (1:1) で3秒				
温度	40 °C				
UV 検出	191/2 nm、リファレンスなし データレート 10 Hz				
MS 検出	ポジティブイオン化モードの SIM				

表 2. MSD スプレーチャンバとシグナルの設定

パラメータ	設定値				
イオン源	Agilent エレクトロスプレー				
ネブライザ圧力	35 psig				
ガス温度	300 °C				
ガス流量	10.0 L/min				
キャピラリ電圧	+5,500 V				
1 秒あたりのターゲットポイント数	3 Hz				
SIM シグナル(m/z)	 63 ジエチルカーボネート (DEC) 77.1 エチルメチルカーボネート (EMC) 89 エチレンカーボネート (EC) 91 ジメチルカーボネート (DMC) 103 プロピレンカーボネート (PC)、ジメチル 2,5- ジオキサへキサンジオエート (DMDOHC) 117 ジエチル 2,5-ジオキサへキサンジオエート (DEDOHC) 147 ジブロピルカーボネート (DPC) 				
フラグメンタ電圧	70 V				

結果と考察

一般的な電解液成分 EC、DMC、DEC、EMC、PC の分離メソッドを開 発して、分解生成物 DMDOHC と DEDOHC の測定も行いました(表 1)。このメソッドは、優れたピーク形状と分解能を実現するために、 InfinityLab Poroshell EC-C18 カラムを使用します。すべての化合物は、 LC/MSD iQ シングル四重極 MSD により検出されましたが、この際、感 度と選択性を向上させるために SIM モードで操作しました。さまざまな 分析対象物の質量電荷比(*m/z*)は、過去の実験で測定されており、そ の際 MSD の設定は Scan モードでした(示していません)。各化合物の 最適な *m/z* を表 2 に示します。

図2は、7種類の分析対象物の定量に使用されたすべてのSIMトレース を重ね表示した、キャリブレーションサンプルの分離結果を示しています。 電解液溶液には使用されていない関連化合物であるDPCを内部標準と して添加しました。図に示したキャリブレーションポイントには、EC、PC、 DMDOHC、DEDOHCが1ppmレベルで含まれていましたが、DMC、 DEC、EMCは1,000ppmで存在していました。内部標準を、すべての サンプルとキャリブラントに1,000ppmでスパイクしました。

すべての分析対象物は、14 分以内に良好に分離されました。SIM トレースは、データ解析メソッドで定義された化合物の予想リテンションタイムのウィンドウ内でのみ表示されます。構造的に類似しているとはいえ、異なる電解液成分が、適用した条件下で大幅に異なるシグナル応答を示したことは注目に値します。EC、PC、分解生成物 DMDOHC と DEDOHC はすべて、DMC、DEC、EMC の約 1,000 倍高い応答を示しました。このため、EC、PC、DMDOHC、DEDOHC のキャリブレーション範囲は 0.01 ~ 10 ppm とし、DMC、DEC、EMC のキャリブレーション範囲は 10 ~ 10,000 ppm としました。直線検量線は、1/x 重み付けと原点除外で作成しました。表3にキャリブレーション結果をまとめ、各化合物の定量下限(LOQ)と検出下限(LOD)を示しています。



図 2. 選択した分析対象物の検出と定量に使用したすべての SIM シグナルのクロマトグラム重ね表示(1,000 ppm/1 ppm キャリブレーションサンプル)

化合物	直線範囲(ppm)	R²	LOQ (ppm)	LOD (ppm)
EC	0.01 ~ 10	0.999	0.147	0.044
PC	0.01 ~ 10	0.998	0.269	0.081
DMC	10~10,000	0.998	52	15
EMC	10~10,000	0.999	94	28
DMDOHC	0.01 ~ 10	0.996	0.011	0.003
DEC	10~10,000	0.999	31	9
DEDOHC	0.01 ~ 10	0.997	0.008	0.002

表 3.7 種類の分析対象物のキャリブレーション結果。LOD と LOQ は、シグナルの 高さ(それぞれ、ノイズレベルの3倍超と10倍超)に基づいて算出しました。

リチウムイオン電池電解液溶液を3回繰り返して分析しました。一定量の純粋なサンプルを純粋なACNで1:10と1:100に希釈して、直接注入に使用しました。100%ACNのような強溶媒中のサンプルは、分離メソッドを低溶媒強度(例えば、2%ACN水溶液)から開始した場合、歪んだピークが生じて、クロマトグラフィーの品質が低下する可能性があります。この問題を解消するため、1260 Infinity II ハイブリッドマルチサンプラを Agilent Feed 注入モードで操作しました。この注入モードでは、溶媒の流れにサンプルを徐々に注入することにより、ピークの歪みやサンプルのブレークスルーなしに、分析対象物をカラムヘッドに濃縮させます。

図 3 に、新しい電解液サンプルの分離結果を示します。主要成分 EC、 DMC、DEC は、1:100 希釈サンプルにおいて高いシグナル応答で検 出されました (図 3A)。DMC と DEC は、それぞれ 32.6 ± 0.6 % と 40.5 ± 0.5 % で定量されました。EC シグナルはキャリブレーション範 囲を上回っていたため、定量できませんでした。そのため、電解液中の EC パーセンテージは、DMC と DEC の合計と 100 % との差として算出 され、26.9 ± 1.1 % でした。PC と EMC はサンプルで検出できません でした。

1:10 希釈では、すでに少量の分解生成物 DMDOHC と DEDOHC が確 認されていました。これは、サンプルを酸素と水のない環境で取り扱う ことができず、分解が生じたためと考えられます。DMDOHC の濃度は LOQ 未満であり、LOQ の半分に相当する 0.007 ppm として報告されて います。DEDOHC は 0.45 ± 0.03 ppm で定量できました。

電解液溶液の経時変化をシミュレートするため、希釈していない一定量 のサンプルを密閉バイアルに入れ、40 ℃ で保管しました。1 週間後、一 定量のサンプルを採取して、ACN で 1:10 と 1:100 に希釈し、内部標準 をスパイクして分析しました。新しいサンプルで検出された分析対象物は すべて、保管しておいたサンプルでも検出されました。DMC と DEC は、 30.7 ± 1.2 % と 39.7 ± 1.5 % という、わずかに低いレベルで定量され ました。これらの数値は、電解液の一部がすでに分解していることを示し ています。EC は再度キャリブレーション限界値を上回りました。DMC と DEC のレベルはすでに低かったため、今回は濃度の推定は行われません でした。不純物である DMDOHC と DEDOHC の量はストレス保管中に 増加しており、それぞれ 12.3 ± 0.04 ppm と 32.6 ± 0.03 ppm と定量 されました。図 4 は、新しいサンプルと、ストレス条件下で 7 日間経過し たサンプルにおける、異なる電解液成分と分解生成物の濃度を示してい ます。



図 3.1:100 (A) と 1:10 (B) に希釈した電解液サンプルの分離結果。クロマトグラムは、すべてのキャリブレーション済み分析対象物の SIM シグナルとピークラベルを示しています。EC はキャリブレーション範囲を上回ったため定量できませんでした。



図 4. ストレス条件下で7日間保管した前後の電解液成分の平均濃度(N = 3)。 EC、DMC、DEC の量はパーセント表示、DMDOHC、DEDOHC は ppm 表示です。

DMDOHC と DEDOHC (m/z 103 と 117) の定量に使用した SIM シ グナルは、異なるリテンションタイムにおいてさらなるピークを示しまし た。これらの分析対象物でモニタリングされた質量電荷比は、インタクト 分子イオンではなく、これらの有機炭酸塩に特異的なフラグメントを示し ています。したがって、ストレスをかけたサンプルに存在する追加のピー クは、DMDOHC および DEDOHC と同じフラグメントを共有している可 能性が高いと考えられます。LC/MS スキャンメソッドによるストレスを かけたサンプルの別の分析では、追加不純物の推定分子イオンが m/z 239、253、281、295 で検出されました。これらの質量電荷比は、LIB 電解液の代表的な分解生成物として報告されているオリゴマー炭酸塩 $C_2H_6(CH_2)_n(CO_3)_3$ (n = 2、3、5、6) の化学式に適合しています。⁴その 生成経路は DMDOHC と DEDOHC と同じであり、LiPF₆の存在下にお いて、EC とそれぞれ 2 当量の DMC または DEC との反応により形成さ れます。DMC と DEC の両方を含む電解液では、EC、DMC、DEC の混 合反応により、追加のオリゴカーボネートが生成されると予想されます。 ただし、適切な参照化合物がないため、想定されるオリゴマーカーボネー トの同定や定量は実施できませんでした。

結論

このアプリケーションノートで紹介しているメソッドを使用して、リチウム イオン電池電解液の主要成分および分解生成物を定量しました。Agilent InfinityLab Poroshell 120 EC-C18 カラムを使用することにより、すべ ての分析対象物を 14 分以内に分離し、優れたピーク形状と分解能を得 ることができました。Agilent 1260 Infinity II ハイブリッドマルチサンプ ラを Feed 注入モードで操作することにより、強溶媒のサンプル(純粋 なアセトニトリル)に起因するピークの歪みが解消されました。Agilent LC/MSD iQ を使用することにより、分析対象物の SIM シグナルの特 異的なモニタリングによる高感度検出が可能になりました。主要成分 の検出下限は 0.04 ~ 28 ppm でしたが、分解生成物 DMDOHC と DEDOHC は 1 桁 ppb の量で検出できましたこのメソッドは、参照化合 物の入手可能性に応じて、さらなる分解生成物を含むように容易に拡張 できます。

参考文献

- Xing, J.; Bliznakov, S.; Bonville, L.; Oljaca, M.; Maric, R. A Review of Nonaqueous Electrolytes, Binders, and Separators for Lithium-Ion Batteries. Electrochem. Energy Rev. 2022, 5, 14.
- Rapid Analysis of Elemental Impurities in Battery Electrolyte by ICP-OES. Agilent Technologies application note, publication number 5994-1937EN, 2020.
- 3. Quick and Easy Material Identification of Salts Used in Lithium-Ion Batteries by FTIR. *Agilent Technologies application note*, publication number 5994-6243EN, **2023**.
- Schultz, C; Vedder, S; Streipert, B; Winter, M; Nowak, S. Quantitative Investigation of the Decomposition of Organic Lithium Ion Battery Electrolytes with LC-MS/MS. RSC Adv. 2017, 7 (45), 27853–27862.

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。

DE40799732

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2024 Printed in Japan, October 15, 2024 5994-6859JAJP

