

Agilent Cary 630 FTIR 分光光度計による 原材料同定アプリケーション

FTIR 用 Agilent MicroLab ソフトウェアを用いて多様なサンプル タイプの分析に対応する高速かつ容易で信頼性の高いワーク フローを実行



著者

Wesam Alwan and Fabian Zieschang Agilent Technologies, Inc.

はじめに

原材料同定は、多くのアプリケーションおよび業界で広く実施されている重要な品質保証および安全管理分 析です。受け取った原材料を同定して確認することで、汚染されたり偽造されたり不正表示された原材料が生 産工程に進むことを防ぐことができます。分析により確実に、最終製品を本来の目的に必要な品質と純度に適 合させることができます。最終製品の同定を確認することは、製品の包装や表示の誤りを回避することと同等 に重要です。したがって、原材料の同定は、世界の多くの薬局方といった業界規格に記載されている要件です。

フーリエ変換赤外分光光度計(FTIR)は、原材料同定の調査で広く使用されており、この目的のために大半の 薬局方で指定されています。FTIRは、化学物質の同定を可能にする、ユーザーが信頼して操作できる高速で 容易で強力な手法です。 適切な分析手法を用いて未知化合物を同定することで、汚染や生産上の問題を調査できます。この分析は、毒物、火薬類、ストリートドラッグ などの危険な物質や違法な物質を同定してファーストレスポンダーおよ び法執行機関に通報するためにも使用されます。

データが有用であることから、さまざまなサンプルに対して原材料同定 ワークフローを熟練していないユーザーがルーチンで実行しています。 このため、分析は高速でシンプルかつ信頼性が高いこと、そして誤った 解釈なしに結果を容易に理解できることが重要です。

Agilent Cary 630 FTIR 分光光度計は、柔軟性があり、高性能なコン パクトベンチトップ機器です(図 1)。モジュール構成の設計のため、ラ ボはさまざまなサンプルタイプおよび FTIR アプリケーションに合わせ て分析システムを構築できます。Cary 630 FTIR 用の相互に交換可能 なサンプルモジュールには、透過、DialPath、TumbliR、ダイアモンド 減衰全反射 (ATR)、ゲルマニウム ATR、ZnSe ATR、ZnSe 多重反射型 ATR、正反射、拡散反射があります。

Cary 630 FTIR 分光光度計は、Agilent MicroLab ソフトウェアを使 用して制御されます。このソフトウェアは画像インタフェースを使用し て、サンプル導入からレポート作成までの分析の各ステップを通じてユー ザーをガイドします(図 2)。どのサンプリングアクセサリがインストール されているかを自動的に検出し、必要な設定を適用して、各サンプルア クセサリ専用の手順画像をロードします。分析結果は色分けされ、ユー ザーはデータを解釈して何らかの適切な処置をとることができます。



図 1. ダイアモンド ATR サンプリングモジュール付き Agilent Cary 630 FTIR を 用いて、異なる粉末の同定を確立します。

この技術概要は、FTIR による原材料同定での異なる手法を紹介します。 アプリケーション例では、さまざまなサンプルタイプに対して、高速で容 易かつ信頼性の高い原材料同定ワークフローを実行するために、Cary 630 FTIR と MicroLab ソフトウェアをどのように使用できるかを示し ます。



図 2. 直感的な Agilent MicroLab ソフトウェアのワークフローにより、Agilent Cary 630 FTIR 分光光度計を用いて簡単に答えにたどり着くことができます。また、画像ガイド 付きソフトウェアは、トレーニングの必要性を低減し、ユーザーのミスのリスクを最小に抑えます。

未知の物質の同定方法およびサンプルの同 定の確認方法

ライブラリの検索

物質にはそれぞれ固有の IR スペクトル (光学異性体および長鎖アルカン 同族体などのいくつかの化合物は除く) があり、物質の指紋と考える ことができます。図 3 に示すように、未知のサンプルの IR スペクトルを 既知化合物のスペクトルライブラリと比較することにより、未知サンプル を同定できます (未知のサンプルがスペクトルライブラリにある場合)。 スペクトルライブラリ内の化合物のスペクトルを用いた「指紋」一致によ る未知の物質の同定は肯定的同定と呼ばれます。

スペクトルライブラリには、数十、数百、時には数千のスペクトルが含ま れている可能性があります。最も良く一致するものを探すために、すべて のスペクトルに対してマニュアルで比較することは多くの場合、困難です。 このため、最も良く一致するものを自動的に同定するための検索アルゴ リズムが開発されています。原材料の同定の確認には、非常に多くの スペクトルを含む大きなライブラリは不要です。多くの場合、目的のアプ リケーションに対して可能性のある材料のスペクトルのみを含むより小 さなライブラリで十分です。ライブラリのサイズに関係なく、ライブラリ 検索のアルゴリズムは、「一致」による同定によって原材料の確認を容易 にします。 ヒットクオリティインデックス (HQI) が各ライブラリのアイテムに対して 自動的に計算されます。HQI 値は、測定したスペクトルとライブラリの スペクトルがどの程度一致するかを示します。HQI は多くの場合、原材 料の同定および確認のワークフローにおいて合格/不合格の基準として 使用されます。

Agilent MicroLab ソフトウェアによるライブラリ検索

Cary 630 FTIR 分光光度計およびアジレントのモバイルおよびハンドヘ ルド FTIR システムに付属する MicroLab ソフトウェアは、メソッドベース の手法を使用します。メソッドをセットアップすると、次のセクションに 記載するように、このソフトウェアによって FTIR システムがターンキー ソリューションに変わり、短時間での決定が可能になります。



図 3. 未知サンプルの FTIR スペクトルを既知化合物のスペクトルと比較して、未 知サンプルを同定できます。このワークフローは「ライブラリ検索」と呼ばれ、 Agilent MicroLab ソフトウェアを使用して自動的に実行できます。

Info Type Instrument Libraries Search Custom Fields Reports	
Search Similarity	
Algorithm: 1650.00 to 1600.00 Derivative Algorithm Gap: 1	Add
Minimum Hit Quality (0-100): 80	
Maximum Hits Displayed: 6	
Quality Critical Threshold 🗹 0.85	
Quality Marginal Threshold 🗵 0.90	Remove
Display Metadata 🗌	Remove All
Auto Residual Search	
Enabled	
Top Hit	
O Defined Spectrum:	

図 4. Agilent MicroLab ソフトウェアで、特定の同定アプリケーションに合わせてライブラリ検索メソッドと検索パラメータをカスタマイズできます。 結果の解釈および報告が容易になるように、結果表示の設定もカスタマイズできます。

メソッドのセットアップ

メソッドのセットアップ時に、ライブラリの検索メソッドを同定アプリ ケーションに合わせてカスタマイズできます。

- 複数のライブラリを同時に検索できます。例えば参照標準のある1
 つのライブラリでサンプルを同定し、規格外サンプルのある別のラ
 イブラリで再発する問題を特定することができます。
- 検索アルゴリズムの幅広い選択肢(ユークリッド、絶対値、微分絶対 値、最小二乗、微分最小二乗、相関、微分相関、類似、微分類似、 拡張相関)を利用できます。
- スペクトル範囲を分析から除外できます。
- 最小 HQI およびヒットの最大数を設定することで、関連するライブ ラリのヒットのみが提供されます。

- HQI を基に色分けされた分析結果を使用して信頼レベルを決定し、結果の解釈を助け、ミスを招く見落としを低減できます。
- その他の化合物固有の情報(危険情報など)を提供できます。

分析の実行

MicroLab は画像を利用して、サンプリング (例えば、固体サンプルに対す る ATR プレスの使用方法) やクリーニング手順など、分析の各ステップで ユーザーをナビゲートします (図 5)。ソフトウェアのホーム画面の**スタート** ボタンをクリックすると、分析が開始されます。

データ取り込み後に、ソフトウェアはライブラリ検索を自動的に実行し、 図6に示すようにベストなライブラリー致のリストを分かりやすい結果 表示フォーマットで提供します。



図 5. Agilent MicroLab ソフトウェアの画像ガイダンスインタフェースは、Agilent Cary 630 FTIR を使用した分析の実行を簡素化し、 トレーニングの必要性を低減します。ソフトウェアは、機器の使用時にオペレータが見ているものと同じ画像を表示します。データ取り 込みの後、図 6 および 10 で示されるように、結果が分かりやすいフォーマットですぐに表示されます。ユーザーが結果を確認した後、 ソフトウェアでは次の分析が開始され、連続的な分析ワークフローが確保されます。



図 6. Agilent MicroLab ソフトウェアは、ライブラリ検索メソッドの結果画面が分かりやすく、経験が浅いユーザーでも確認して結果を解釈できるようになっています。左:ライブ ラリヒットデータが、リストのみで、またはリストとスペクトルで表示されます。右:危険情報やファーストレスポンス情報を含む化学物質メタデータのオプション画面も利用できます。

ライブラリとライブラリ管理

ライブラリ検索の品質は、ライブラリの品質に大いに依存します。アジレントは、すぐに使用できる、MicroLab ソフトウェアと使用できるアプリケーション固有のライブラリを豊富に提供しています。

また、MicroLab ソフトウェアで、スペクトルライブラリを容易に作成、 保守、管理することもできます(図 7)。

新しいライブラリを数秒で作成することができます。作成時にも、その他のいつでも、結果画面からすぐにスペクトルをライブラリに追加できます。

	User: Method:	User New method - ol	ld design	
LibraryNar Agilent Ela Agilent Po API and ex	me astomer Oring and S Ilymer Handheld ATR xcipient.lib	eal Handheld ATR Library L Library.lib	Contaif /.lib	Create Library X
Library Cor	mment:			Comment This library includes new compounds
Index 0 1 2	Deleted?	CAS# 10016-20-3 15687-27-1 63-42-3	Entry Nar Cyclodext Ibuprofen Lactose -	Library Spectra are: Baseline Corrected ATR Spectra Y Type: Lib Bytes: Absorbance V 2 Bytes V
	3500			Auto-detect from first library entry
	_	Home Add To) ,	Create Cancel

図7. Agilent MicroLab ソフトウェアではライブラリ管理は高速かつ簡単です。新規ライブラリは数秒で作成でき、必要に応じて ライブラリの内容を更新できます。

アプリケーション例

アプリケーション例1:医薬品の原材料の同定確認

ダイアモンド ATR サンプリングモジュール付き Cary 630 FTIR を使用し て、医薬品有効成分 (API) でよく使用される 3 成分、サリチル酸、イブ プロフェン、ニコチン酸および、賦形剤 (α-シクロデキストリン) 参照物 質の IR スペクトルを収集しました。MicroLab ソフトウェアでスペクトル ライブラリが作成され、収集した参照スペクトルがこのライブラリに追加 されました。このスペクトルライブラリをライブラリ検索メソッドで使用し、 図 8 で示すパラメータを使用しました。 次に、この原材料同定確認のメソッドを使用して、サリチル酸、ニコチン酸、 イブプロフェンおよび α-シクロデキストリンの異なるバッチを分析しま した。MicroLab ソフトウェアで、すべてのテストされた原材料の同定が 正しく確認されました。HQI はそれぞれ、0.99238 (サリチル酸)、0.99775 (ニコチン酸)、0.98641 (イブプロフェン)、0.98675 (α-シクロデキスト リン) でした。ここで、1 は、類似性検索アルゴリズムを使用した最も高 い理論値です (図 9)。

Info Type Instrument Libraries Search Custom Fields Reports	
Search Similarity Exclude Search Regions (cm-1):	
Derivative Algorithm Gap: 1	Add
Minimum Hit Quality (0-100): 50	
Maximum Hits Displayed: 6	
Quality Critical Threshold 🗹 0.90	
Quality Marginal Threshold 🛛 0.95	Remove
Display Metadata	Remove All
Auto Residual Search	
Enabled	
Top Hit	
O Defined Spectrum:	

図 8. 医薬品の API の同定を確認するために使用したライブラリ検索パラメータ。品質重要しきい値と 品質限界しきい値の設定により、結果表示での色分けが決定されます。



図 9. Agilent MicroLab ソフトウェアではライブラリ検索結果で色分けが適用されます。色分けにより、 結果の解釈が簡素化され、ミスを招く見落としが低減されます。スクリーンショットでは、α-シクロデキス トリンの分析結果が緑色で表示され、ライブラリスペクトルと適切に一致することを示しています。

アプリケーション例2:ポリマーサンプルの同定

ダイアモンド ATR サンプリングモジュール付き Cary 630 FTIR を使用して、ポリプロピレン、ポリカーボネート、塩素化ポリエチレン、ポリエチレン テレフタレート、ポリ塩化ビニルを含む 5 つの異なるプラスチック材料を 分析しました。

Agilent FTIR スペクトルライブラリATR Polymers and Polymer Additives (部品番号 G8045AA、オプション 106) を Agilent MicroLab ソフトウェアで選択しました。このスペクトルライブラリには、代表的なポ リマー、プラスチック、ポリマー添加剤、可塑剤、充填剤のスペクトルが 7,974 個含まれています。

スペクトルライブラリを検索するために類似性アルゴリズムを使用し、 HQI 結果によって示されるように、5つのポリマーサンプルすべてが適切に 同定されました。HQI の値は、0.98315 (ポリプロピレン)、0.97298 (ポリ カーボネート)、0.97212 (塩素化ポリエチレン)、0.96238 (ポリエチレン テレフタレート)、および 0.98360 (ポリ塩化ビニル) でした。ここで、1 は、 類似性検索アルゴリズムを使用した最も高い理論値です(図 10)。



図 10. ダイアモンド ATR サンプルモジュール付き Agilent Cary 630 FTIR を使用 して得られる色分けされた一致結果が短時間で明確に表示されます。ルーチン、定性、 材料同定メソッドにより、高い信頼性でポリマーの種類を自動的に同定できます。

スペクトルのマニュアルでの比較

既知の原材料の同定は、IR スペクトルのマニュアル比較によっても確認 することができます。サンプルのスペクトルは参照物質のスペクトルと目視 で比較するか、または測定されたピークがラベル付けされこれを文献値 と比較します。

通常、ほとんどのスペクトル情報は、IR スペクトルの指紋領域(約400~ 1,500 cm⁻¹)内にあります。MicroLab ソフトウェアを使用すると、スペ クトルのエネルギーがより高い領域の波数スケールを圧縮しつつ、指紋 領域の波数スケールを拡張できます(図 11)。スペクトルを拡張すること で、より関連する指紋領域をより詳しくかつより容易に調査できます。



図 11. 物質の IR スペクトルは、官能基の存在など原材料の組成に関して価値ある 情報を提供します。多くの場合、大半の関連するバンドは、IR スペクトルの指紋領域 (約 400 ~ 1,500 cm⁻¹)内で見つけられます。Agilent MicroLab ソフトウェアで は、波数のスケーリング係数を適用でき、対象のスペクトル範囲のより詳細な分析 を可能にします。

アプリケーション例 3:マニュアルでのスペクトル比較による API の原材料同定の確認

透過サンプリングモジュール付き Cary 630 FTIR を使用し、リファンピン (抗生物質 API) を含む臭化カリウム (KBr) ペレットの IR スペクトルを 収集しました。4,000 ~ 2,000 cm⁻¹ の範囲にはスペクトル情報がほとん ど観察されず、2,000 ~ 600 cm⁻¹ の領域に大部分のバンドが存在しま した。MicroLab ソフトウェアのグラフィック結果画面で、より重要な指紋 領域を (係数を2として) 拡張し、より高い波数のスペクトルは圧縮しま した (図 12)。 スペクトルを拡張することで、サンプルスペクトルを国家薬典委員会が 発行する Atlas of infrared spectra of drugs に記載されている参照 スペクトルと容易に比較できました。関連するバンドにはラベルが付けられ、 スペクトル、およびすべてのラベル付きピークの表を含む PDF レポート がソフトウェアで自動的に生成されました(図 13)。



図 12. リファンピンの IR スペクトル。スペクトルを確認できるように、波数がより高い領域と比べて係数2で指紋領域を拡張し、関連するバンドにラベルを付けました。



図 13. Agilent MicroLab ソフトウェアの結果ウィンドウからすぐに PDF レポートを生成できます。レポートはフルカスタマイズ可能で、すべての関連する情報、例えば、サンプル名やオペレータ ID および、拡大されたスペクトル、フルスペクトル、ピークリストを含む分析の詳細を含めることができます。

結論

FTIR による原材料の同定は多くの業界で重要な分析です。この技術概要で は、アプリケーションに適したサンプリングモジュールとカスタマイズされた メソッドを使用し、多様なサンプルタイプの原材料の同定において Agilent Cary 630 FTIR に柔軟性があることが示されました。ワークフローを通じて 段階的にガイダンスを提供する Agilent MicroLab ソフトウェアを使用し、 メソッド開発を簡素化しました。アプリケーション例では、ライブラリ検索や スペクトルのマニュアル比較による原材料の同定が、ソフトウェアによってど のように容易になるかを示しました。

直感的な、画像ガイド付き MicroLab ソフトウェアは短時間で簡単に習得で きるため、トレーニングの必要性およびオペレータのミスが低減し、レポート 作成時間が短縮するため、原材料の品質管理試験の生産性が向上します。

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。

DE22553965

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2022 Printed in Japan, July 20, 2022 5994-4992JAJP

