

GC/Q-TOF と Low-Energy El による 結核感染マウス肺組織の代謝変化の分析



著者

M^a Fernanda Rey-Stolle¹, Vineel P. Reddy², Santiago Angulo¹, Adrie J.C. Steyn^{2,3,4}, Sofia Nieto⁵, Nathan Eno⁵, and Coral Barbas¹

¹ CEMBIO, Facultad de Farmacia, Universidad CEU San Pablo Madrid, Spain

² Department of Microbiology, University of Alabama at Birmingham Birmingham, AL

- ³ KwaZulu-Natal Research Institute for TB and HIV (KRITH) Durban, South Africa
- ⁴ UAB Center for Free Radical Biology, University of Alabama at Birmingham Birmingham, AL
- ⁵ Agilent Technologies, Inc. Santa Clara, CA

概要

結核 (TB) の世界的蔓延による罹患者の増加は甚大で、2014 年だけでも新たな TB 発症者が 960 万 人、この疾患による死亡者は 150 万人と見積もられています [1]。メタボロミクスを用いて TB バイオ マーカーを同定することにより、この疾患に対する理解を深めることができます。この研究では結核 菌 (Mtb) 感染のマウスモデルを用いて、非感染および感染肺組織の代謝プロファイルを特定しました。

感染に関係する新しい病態生理学的経路と TB のバイオマーカーを同定するために、非感染肺組織 と感染後 9 週間が経過した肺組織を用いて、ノンターゲットメタボロミクス分析を実施しました。最 初の化合物アノテーション後、Low-Energy El データを使用して分子イオンを確認し、暫定的に同定し た化合物と未知化合物それぞれの分子式を同定しました。



実験方法

気管内経路を通して 5×10^4 CFU の結核菌 (Mtb) H37Rv を投与し、マウス を感染させました。0-Xトキシ化した後にトリメチルシリル化し、肺組織 の乾燥抽出物を誘導体化しました。Agilent 7890B GC システム、高分解能 Agilent 7250 GC/0-TOF、および Low-Energy イオン化が可能な El イオン源を 組み合わせて、GC/MS 分析を実施しました (図 1)。Agilent 7250 システム に Low-Energy El イオン源を追加することにより、高分解能 (m/2 272 にお いて 25,000) が達成されて質量精度が向上し、ダイナミックレンジが広が ります。表1に、分析条件を示します。

最初の化合物同定で Fiehn.L RI ライブラリを用いる際の化合物同定を容易にするために、リテンションタイムロッキング (RTL) Fiehn メソッドを使用 しました。さらに別の化合物を同定するために、米国国立標準技術研究 所のライブラリ (NIST.L) と精密質量メタボロミクス用 PCDL を使用しまし た。Agilent MassHunter Unknowns Analysis B.08.00 の SureMass 機能を用い て、フィーチャー検出を実施しました。Agilent Mass Profiler Professional (MPP) バージョン 13.0 で統計解析を実施しました。最後に、MPP の拡張ツール である Pathway Architect を用いて、TB 感染に関係する生化学的経路を同 定しました。



🗵 1. Agilent 7250 GC/Q-TOF

表 1. Agilent 7250 GC/Q-TOF 分析条件

パラメ-タ	設定
カラム	Agilent DB-5MS、 30 m × 0.25 mm、0.25 μm、 DuraGuard、10 m
注入量	1 μL
スプリット比	10:1
スプリット/スプリットレス注入口 温度	250 ° C
オーブン温度プログラム	60°Cで1分、 10°C/minで325°Cまで上昇、 9.5分維持
キャリアガス	ヘリウム、1 mL/min 定流量
トランスファライン温度	290 ° C
イオン化モード	標準 El、70 eV Low-Energy El、17 eV、15 eV、および 12 eV
イオン源温度	200 ° C
四重極温度	150 ° C
質量範囲	$50\sim950~m/z$
スペクトル採取レート	5 Hz

結果と考察

実験の設定とフィーチャー検出

感染に関係する新しい病態生理学的経路と TB のバイオマーカーを同定 するために、非感染肺組織抽出物と感染後 9 週間が経過した肺組織抽 出物を用いて、ノンターゲットメタボロミクス分析を実施しました。

Unknowns Analysis で実施したフィーチャー検出とライブラリ検索 (図 2) の 結果を、さらに MPP で処理するために .CEF ファイルとしてエクスポートし ました。

主成分分析によるクラスタリング

MPP での主成分分析 (PCA) により、データのクラスタリングを評価しました。非感染対照群と感染組織の間で、完全に異なる特徴的なクラスタが 形成されました (図 3)。



図 2. Agilent MassHunter Unknowns Analysis (例として PCDL を使用) で実施したフィーチャー検出とライブラリ検索



図 3. PCA プロット。非感染対照群 (UC、青色の円) と感染後 9 週間が経過した (9W、赤色の円) 肺組織の間で観察された特徴的なクラスタ

差分解析

感染マウスと非感染マウス間の肺組織メタボロームにおける有意な変 化を MPP で詳細に評価するために、倍率変化解析 (図 4) とヒートマップ (図5)を使用しました。多数の代謝物、特にアミノ酸と核酸塩基のプロファ イル内で変化が観察されました。さらに、イタコン酸とキヌレニンのプロ ファイル内でも変化が検出されました。

.... -1 5 6 イタコン酸 . . in the 5 -log₁₀(補正後の p-値) 4 3 2 .1 all a 1 11 120 12 嚎 CONTRACTOR OF OGER ROOM COLUMN TWO IS NOT 0 1000 (CHERNER) 0 10 -20 20 -10 0 log₂(倍率変化) Select pair [UC] Vs [9W]

Control Group [9W] 2.0





図 5. 非感染肺組織と感染肺組織間の規制対象代謝物の差を示したヒートマップ

TB 感染に関係している可能性のある生化学的経路の同定に対しても Pathway Architect を用いました。プリンとピリミジン代謝、さらに NAD 生 合成 || の経路は最も重要なものの1つであり、図6にその例を示してい ます。



図 6. 経路解析の結果例: NAD 生合成 II

未知化合物の同定と暫定的に同定した化合物の確認

最初の化合物アノテーションおよび MPP での差分解析後、Low-Energy El スペクトルを使用して分子イオンを確認し、暫定的に同定した差分化合物と未知化合物の分子式のそれぞれを同定しました (図 7)。

未知化合物の構造を解析するための最初のステップでは、Low-Energy El により分子イオン (M^+) を同定しました。さらに、 M^+ の絶対アバンダンス をプリカーサとして最大限に使用するために、最適なイオン化エネルギー において MS/MS を取得しました。その後、Qual で Find by Targeted MS/MS アルゴリズムを用いて、未知化合物の MS/MS スペクトルを抽出しました。 その結果を、Molecular Structure Correlator (MSC) で評価しました (図 8)。



図 7. Low-Energy El による分子イオン (M+) の確認



60 70 80 90 100 110 120 130 140 150 160 170 180 190 200 210 220 230 240 250 260 270 280 290 質量電荷比 (m/z)



図 8. Low-Energy El、MS/MS、および Molecular Structure Correlator (MSC) による未知化合物の構造解析。 未知化合物に関して可能性のある構造を図示

結論

今回のノンターゲットメタボロミクス分析では、アミノ酸のプロファイルおよびキヌレニンとイタコン酸のプロファイル内で変化が検出されました。中でも興味深いのは、イタコン酸は一般的に哺乳類の代謝物としては分類されていないにもかかわらず、最近になって、マクロファージベースの免疫応答において一定の役割を果たしているということが明らかになったことです [2]。

参考文献

- World Health Organization (http://www.who.int/mediacentre/factsheets/ fs104/en/)
- 2. Cheryl L. Strelko, et al. J. Am. Chem. Soc. 133(41), 16386-16389 (2011).

詳細

本文書のデータは代表的な結果を記載したものです。 アジレント製品とサービスの詳細については、アジレントのウェブサイト www.agilent.com/chem/jp をご覧ください。 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. 2017 Printed in Japan, July 6, 2017 5991-8199JAJP

