

The Agilent Metabolomics Dynamic MRM Database and Method

Technical Overview

Author

Mark Sartain
Agilent Technologies, Inc.
Santa Clara, California, USA

Introduction

A major challenge in metabolomics is achieving reproducible, robust chromatographic resolution with a single analytical LC/MS method for endogenous cellular metabolites due to their diversity of chemical and physical properties. A more manageable chromatographic solution is to develop pathway-targeted analytical methods for smaller sets of metabolites. Building an optimized targeted LC/MS/MS method is a painstaking task requiring significant time and money to characterize a large number of chemical standards.

Agilent has developed a LC/MS/MS solution for >215 central carbon metabolites that provides robust, reproducible separation of acidic metabolites. The solution includes a curated database with retention times, optimized MS/MS acquisition parameters, and a data acquisition and analysis method. The Metabolomics dMRM Database and Method leverages the high sensitivity and wide dynamic range of the Agilent 6470 Triple Quadrupole LC/MS System, and the proven ionization enhancement capabilities of the Agilent Jet Stream Ionization Source [1].

This application note illustrates the advantages of the Agilent Metabolomics dMRM Database and Method for enabling researchers to easily implement an already-optimized and straightforward targeted analytical method for central carbon metabolites.



Agilent Technologies

Solution Overview

Many chromatographic approaches have been applied to metabolomics analysis, including reversed-phase, normal phase, and HILIC, but all are generally unsatisfactory due to poor selectivity, reproducibility, or retention of polar metabolites. To address these challenges, a highly reproducible and robust ion-pair reversed-phase (IP-RP) chromatographic method has been developed to provide separation of anionic and hydrophobic metabolites. The chromatographic method leverages the robustness of Agilent ZORBAX Extend C18 columns with the ion-pairing agent tributylamine (TBA). Normally a nonvolatile quaternary amine would be used for ion pair chromatography, however, quaternary amines interfere with electrospray ionization. TBA is a volatile tertiary amine amenable to electrospray ionization that functions as an ion pair reagent and facilitates reproducible retention of acidic metabolites. Nonacidic metabolites still maintain conventional reversed-phase selectivity in the presence of TBA. The ion-pairing LC/MS method described here enables simultaneous analysis of many metabolite functional classes, including amino acids, citric acid cycle intermediates and other carboxylic acids, nucleobases, nucleosides, phosphosugars, and fatty acids.

Instrumentation

Coupling the Agilent 1290 Infinity Series UHPLC to an Agilent 6470 Triple Quadrupole LC/MS creates a sensitive and highly selective system for the detection of targeted metabolites (Figure 1). To achieve high sensitivity and selectivity, a triple quadrupole is operated in multiple reaction monitoring (MRM) mode. MRM mode specifies for each compound a predetermined precursor ion and collision-induced product ion pair, which is termed a transition. In complex samples, a second transition is highly desirable, if possible, to increase confidence in compound identification. Including chromatographic retention time of the transitions provides orthogonal confirmation of compound identification.

When analyzing many compounds in a single LC/MS analytical method, the number of concurrent MRM transitions increases, which reduces measurement (dwell) times and impacts limits of detection (LODs). With retention time information for each compound, MRM transition lists can be dynamically created throughout an LC/MS run using a window around the expected retention times. In this way, compounds are only monitored while they are eluting from the LC, thus improving LODs and permitting more metabolites to be measured. In Agilent MassHunter Acquisition Software, using retention time windows is called dynamic MRM (dMRM) mode [2]. This solution uses a dMRM method to maximize LODs for central carbon metabolites.



Figure 1. Agilent 1290 Infinity II UHPLC system with an Agilent 6470 Triple Quadrupole Mass Spectrometer.

This analytical method is intended for use with complex samples where matrix components may impact method performance. The LC/MS configuration for this solution includes a guard column and a switching valve configuration that permits backflushing of both the column and guard column following each analytical run. This very effectively clears potential matrix contaminants from the guard column, and provides consistent high-quality peak shapes and superior long-term retention time stability.

Methodology

The IP-RP chromatography provides effective separation of anionic and hydrophobic metabolites. This method uses a gradient between solvent A (97:3 water/methanol with 10 mM tributylamine + 15 mM acetic acid) and solvent B (methanol with 10 mM tributylamine + 15 mM acetic acid). We choose the Agilent ZORBAX Extend C18 column as it is very stable at the pH used with this method. More detailed information about the chromatography and solvent preparation can be found in the Metabolomics Dynamic MRM Method Setup Guide (G6412-90002).

The Agilent advanced triple quadrupole Autotune program optimizes all noncompound-dependent instrument settings. Due to the different physicochemical properties of metabolites, there are specific mass spectrometer parameters that must be optimized for each compound to provide the best performance.

On the Agilent system, this includes fragmentor voltage, which optimizes transmission of the precursor ion into the mass spectrometer, and collision energy, which provides maximum intensity for a specific product ion. The Agilent Metabolomics dMRM Database provides optimized parameters for all metabolites. The choice of primary transitions, retention times, collision energies, and fragmentor voltages included in the database were developed in collaboration with the laboratory of Adam Rosebrock (University of Toronto), a leading group in the metabolomics field. These optimized parameters were the result of extensive analyses of chemical standards to generate results of the highest quality.

The database links to both MassHunter Acquisition Software and MassHunter Optimizer [3]. MassHunter Optimizer can be used to add additional metabolites to the database if needed. Users can select compounds from the database to rapidly create a customized analytical method for their specific needs. Figure 2 shows the MS QQQ Acquisition setup tab for controlling the Agilent 6400 Series Triple Quadrupole LC/MS system. The database contains metabolite information such as compound name, CAS number, and the optimized acquisition parameters for each transition. In addition, retention times with retention time windows are included for Dynamic MRM mode, which are valid if the user is implementing the same hardware configuration and LC method supplied with the database. Using the Agilent Database Browser (Figure 3), compounds with their optimized acquisition parameters can be selected for import into the acquisition method.

The screenshot displays the Agilent MassHunter Acquisition software interface, specifically the QQQ tab. The interface is divided into several sections:

- Properties:** Includes fields for Tune file (tunes.TUNE.XML), Stop time (2.4 min), Ion source (AJS ESI), and Time filtering (Peak width 0.04 min).
- Time segments:** A table with columns for #, Start Time, Scan Type, Div Valve, Delta EMV (+), Delta EMV (-), and Stored. Row 1 shows a Dynamic MRM scan at 0 minutes.
- Acquisition:** A table with columns for Compound Group, Compound Name, MS2 Res, and Ret Time (min). A context menu is open over the table, with 'Import from Database Browser...' highlighted. Other menu items include Insert Row, Append Row, Delete Row, Sort, Update DMRM Method..., Edit DMRM Method..., Calibrate DMRM Method..., Cut, Copy, Paste, Paste from Clipboard, Fill Down, and Fill Column.
- Dynamic MRM Parameters:** Shows Cycle Time set to 500 ms.

Figure 2. Agilent MassHunter Acquisition QQQ tab showing the **Import from Database** function.

Database Browser

File Edit View

Search/Filter Import List

Show All Records Search/Filter

Filter Compounds

Enable Filters

Optimized Compounds

Date From 11/19/2015 To 06/04/2010

Group Name Project Name

Polarity Positive Model

Method

Search Compounds

Search Text

Select Columns

- Project Name
- Compound Name
- Formula
- MW
- Groups
- CAS
- Chemical Classes

Match entire word for each string

Select Transitions

Select top 1 ranked transitions

Primary transitions

Secondary transitions

Select Transitions

Set primary and trigger flags

Set top 2 ranked transitions as primary

Set Primaries and Trigger

Rank transitions by

Abundance

Response Factor

<input type="checkbox"/>	Compound Name	Formula	Precursor	Product	Frag	CE	CAV	RT	RT Window
<input checked="" type="checkbox"/>	L-Methionine	C9H11NO2S	148	100	70	4	4	1.667	1
<input checked="" type="checkbox"/>	L-Methionine	C9H11NO2S	148	47.2	70	8	4	1.667	1
<input checked="" type="checkbox"/>	L-Histidine	C6H9N3O2	154.1	137	102	12	4	1.087	1
<input checked="" type="checkbox"/>	L-Histidine	C6H9N3O2	154.1	93.1	102	16	4	1.087	1
<input checked="" type="checkbox"/>	L-Phenylalanine	C9H11NO2	164.1	147	100	8	4	3.9	1
<input checked="" type="checkbox"/>	L-Phenylalanine	C9H11NO2	164.1	103	100	16	4	3.9	1
<input checked="" type="checkbox"/>	L-Glutamine	C9H17NO3	145.1	127	94	8	4	1.242	1
<input checked="" type="checkbox"/>	L-Glutamine	C9H17NO3	145.1	109	94	8	4	1.242	1
<input checked="" type="checkbox"/>	L-Tryptophan	C11H12N2O2	203.1	116	116	16	4	6.667	1
<input checked="" type="checkbox"/>	L-Tryptophan	C11H12N2O2	203.1	74.1	116	12	4	6.667	1
<input checked="" type="checkbox"/>	L-Threonine	C4H9NO3	118	74.1	72	8	4	1.25	1
<input checked="" type="checkbox"/>	L-Serine	C3H7NO3	104	74.1	70	8	4	1.233	1

Current Database: D:\MassHunter\Databases\QQQ_dMRM_Metabolomic_DB

Add to Import List Import Close

Figure 3. Database Browser view of Metabolomics dMRM Database.

Key to the LC/MS analytical method described here are the compound retention times in the database. The retention times provide added confidence in assigning the MRM signals to the correct metabolite in a complex biological matrix, and enable Dynamic MRM Acquisition. The retention time consistency of the IP-RP method allows narrow retention time windows, enabling the analysis of larger numbers of metabolites with better LODs.

A default acquisition method (.m) file is provided with the database for seamless implementation of the full Dynamic MRM method targeting all metabolites. The method file provides all of the required instrument parameters, many of which are essential to replicate the retention times specified in the database. The Agilent Jet Stream source settings provided in the analytical method were optimized for efficient ionization for the broad range of metabolites. The cycle time was chosen to maximize sensitivity and ensure sufficient data points for reproducible chromatographic peak integration. Figure 4 shows the overlapped MRM chromatograms of a metabolite standard mixture acquired with the default acquisition method.

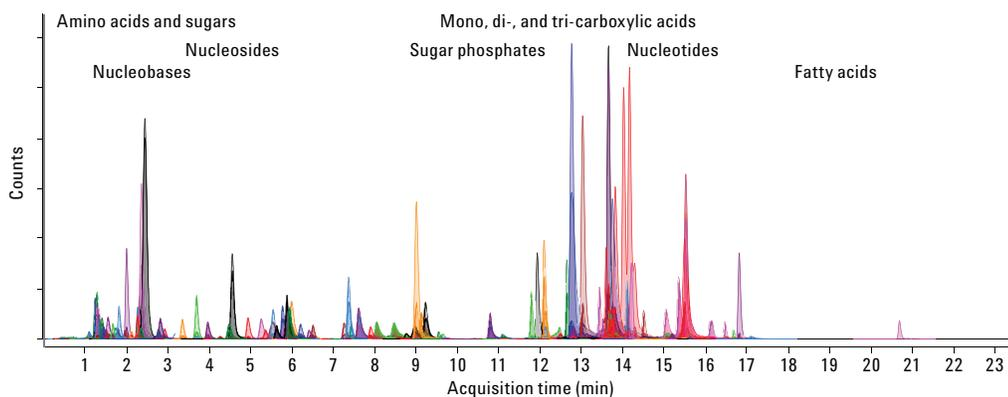


Figure 4. Overlaid MRM chromatograms of more than 100 metabolite standards at 5 ng on-column.

Data Analysis Workflow

The Metabolomics Dynamic MRM (dMRM) Database and Method includes a MassHunter Quantitative Analysis method for processing the LC/MS data. For large, multicomponent batches, MassHunter Quantitative Analysis Software facilitates batch review of the data by including helpful features such as compound-at-a-glance data review (Figure 5), dynamically linked results, and parameterless integration algorithms.

The analytical method described here has been designed to enable relative comparison studies. The data analysis method could be quantitative if metabolite calibration curves were added by the user. In addition, the method, database, and workflow could be extended by the user to include spike-in controls and transitions for isotopically-labeled internal standards to achieve better quantitation.

Results from the Metabolomics Dynamic MRM Database and Method are supported in Agilent comprehensive metabolomics workflows. To achieve this, a report is automatically generated (Figure 6) from MassHunter Quantitative Analysis Software for import into the Agilent multivariate analysis software, Mass Profiler Professional [4]. The metabolite names, integrated peak abundances, and the chemical identifiers (CAS) are included in the import for multivariate sample comparisons as well as subsequent pathway analysis. The differential results can be used in the Pathway Architect module of Mass Profiler Professional to identify affected pathways for biological interpretation (Figure 7). Pathway Architect supports three major sources of pathway content: WikiPathways, BioCyc/MetaCyc, and KEGG [5]. For increased biological understanding, the targeted metabolomics results can be combined with other omics data (proteomics, transcriptomics) for a multi-omics analysis within Pathway Architect.

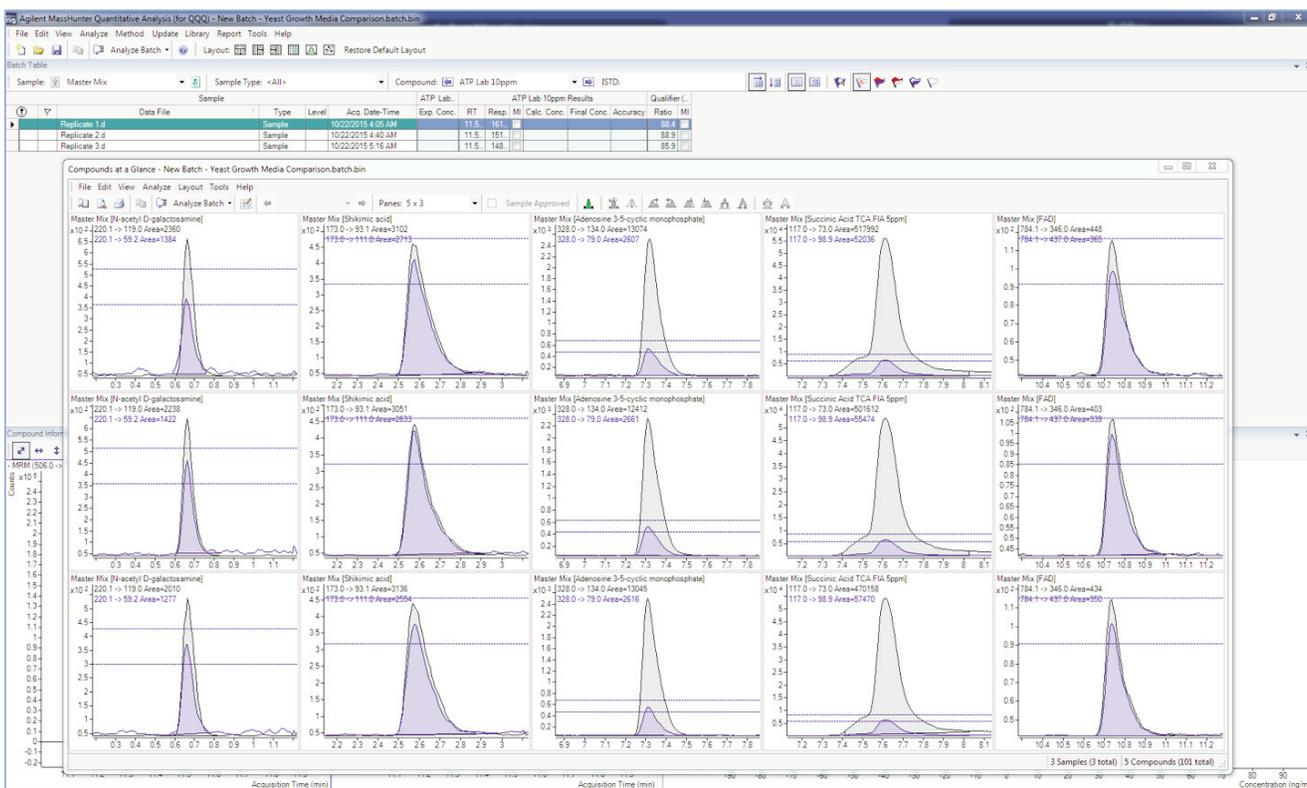


Figure 5. Compounds-at-a-glance view within Agilent MassHunter Quantitative Analysis Software, displaying five compounds (columns) across three samples (rows).

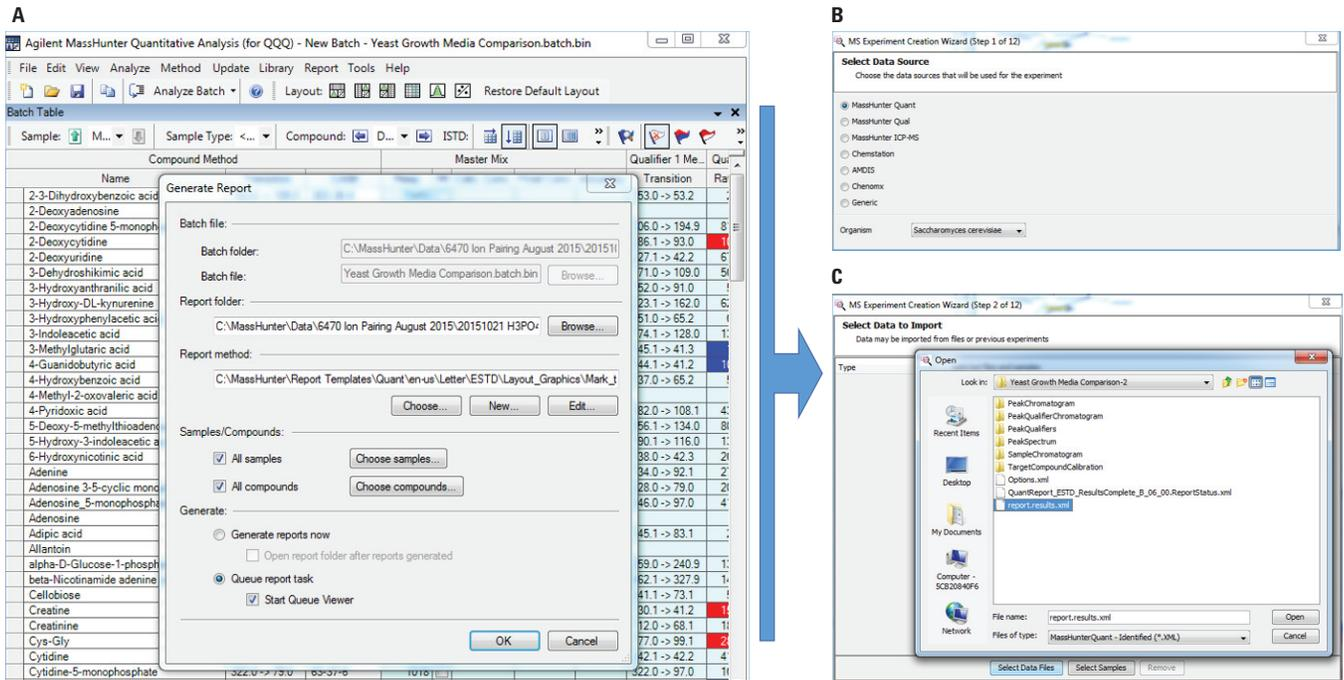


Figure 6. Agilent MassHunter Quantitative Analysis Report generation and import into Agilent Mass Profiler Professional. A report is easily generated from the results of a large sample batch (A), and the results are imported into Mass Profiler Professional after the data source type is selected (B) and the location of the Quant report file is specified (C).

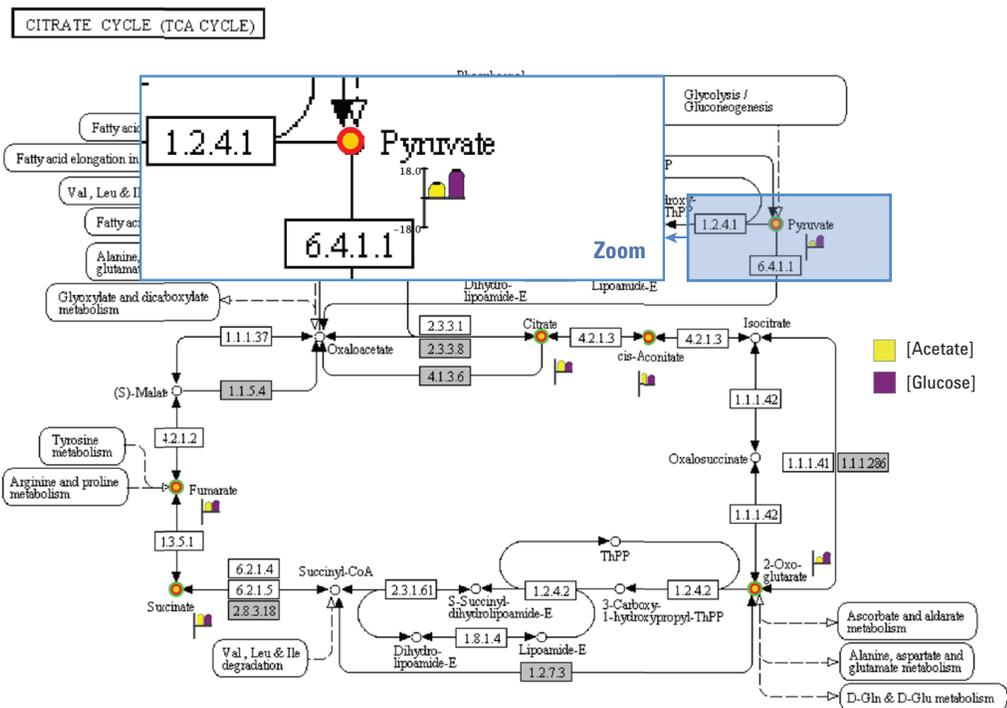


Figure 7. The Pathway Architect module of Agilent Mass Profiler Professional, showing the Yeast TCA pathway from KEGG. Metabolites matching experimental data are highlighted. Data is visualized on these pathways with heat strips indicating the normalized relative abundances between sample groups, in this case yeast grown cultivated on acetate or glucose.

Conclusions

The Agilent Metabolomics Dynamic MRM (dMRM) Database and Method was developed for targeted LC/MS analysis of central carbon metabolites in complex biological matrices. The LC/MS method is optimized for an Agilent 1290 Infinity II UHPLC with an Agilent 6470 Triple Quadrupole Mass Spectrometer. The combined UHPLC-MS/MS method and acquisition database provides; robust ion-pairing reversed-phase chromatography optimized for challenging acidic metabolites, a curated database including metabolite-specific LC/MS/MS transitions and retention times, and a default data analysis method, which is supported in Agilent comprehensive metabolomics workflows.

The Metabolomics Dynamic MRM (dMRM) Database and Method provides rapid setup and excellent results eliminating the need for analytical method development.

References

1. S. Buckenmaier, *et al.* "Instrument Contributions to Resolution and Sensitivity In Ultra High Performance Liquid Chromatography Using Small Bore Columns: Comparison Of Diode Array and Triple Quadrupole Mass Spectrometry Detection" *J. Chromatogr. A* **1377**, 64-74, (2015).
2. P. Stone, *et al.* New Dynamic MRM Mode Improves Data Quality and Triple Quad Quantification in Complex Analyses. *Agilent Technologies Technical Overview*, publication number 5990-3595 (2009).
3. MassHunter Optimizer Software for Automated MRM Method Development Using the Agilent 6400 Series Triple Quadrupole Mass Spectrometers. *Agilent Technologies Technical Overview*, publication number 5990-5011 (2010).
4. Agilent Mass Profiler Professional Software: Discover the Difference in your Data. *Agilent Technologies Brochure*, publication number 5990-4164 (2014).
5. Agilent Pathway Architect: From Discovery to Insight. *Agilent Technologies Brochure*, publication number 5991-1459 (2014).

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

For Research Use Only. Not for use in diagnostic procedures.

This information is subject to change without notice.

© Agilent Technologies, Inc., 2016
Printed in the USA
March 10, 2016
5991-6482EN



Agilent Technologies