

Use of Charged Aerosol Detection as an Orthogonal Quantification Technique for Drug Metabolites in Safety Testing (MIST)

Gregory Malek,¹ Chris Crafts,² Marc Plante,² Matthew Neely,¹ and Bruce Bailey²

¹Thermo Fisher Scientific, Bannockburn, IL, USA; ²Thermo Fisher Scientific, Chelmsford, MA, USA



Abstract

The U.S. FDA guidelines on drug metabolites in safety testing (MIST) published in 2008 requires that the relative quantitation of human metabolites be obtained as soon as feasible in the drug development process. The goal is to ensure that at least one of the animal models used is not only producing the same metabolites found in humans, but at the same relative abundance. Identification of drug metabolites is typically accomplished using various LC-MS approaches. However, metabolite quantification is often difficult to accomplish with MS techniques alone. In some cases, the addition of UV detection is sufficient for metabolite quantification. But this requires that both parent drug and its metabolites possess similar and sufficiently active chromophores, which is not always the case. Charged aerosol detection can detect any nonvolatile compound, typically with low ng sensitivity and similar response independent of chemical structure.

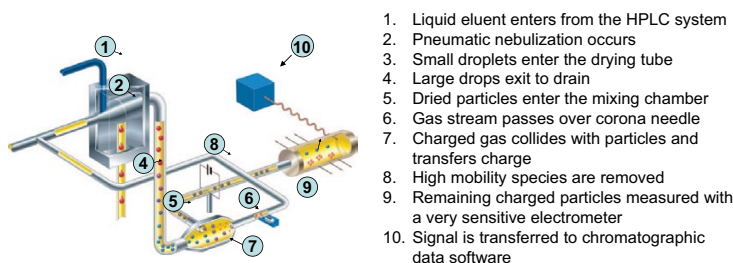
Presented here are the results from two test cases used to evaluate the usefulness of charged aerosol detection for MIST: buspirone (which has a strong UV chromophore) and erythromycin (which has a weak UV chromophore). These drugs were analyzed directly following incubation with human liver microsomes (drugs at $\leq 100 \mu\text{M}$). The five major metabolites for buspirone and the four major metabolites for erythromycin were identified using an a Thermo Scientific LTQ Orbitrap mass spectrometer and quantified by charged aerosol detection. The range, linearity, and sensitivity of this approach are discussed. Charged aerosol detection provides a complementary technique to those already being employed in industries capable of extending the range of in vitro drug metabolites that can be monitored during the drug development stage.

Introduction

Interest in metabolite and trace impurity analysis by the pharmaceutical industry is intensifying due to concerns with mass balance studies, regulatory commitments in reporting active pharmaceutical ingredient (API) impurities, MIST, and cleaning validation of manufacturing equipment. Most often an analytical requirement for accurately reporting the level of metabolites or impurities is to obtain reference standards or use a radiolabeled drug approach. Because many of these standards are unavailable and the radiochemical approaches are time consuming and expensive, quantification of drug metabolites can often be difficult in early development stages. The situation is further exacerbated because several types of HPLC detectors, such as UV or evaporative light scattering detection (ELSD), either do not provide uniform response across the target analytes or lack the sensitivity to detect these compounds. Although electrospray ionization (ESI) MS techniques are extremely sensitive and powerful approaches, ionization efficiency can vary between metabolites and lead to quantification issues.

The Thermo Scientific Dionex Corona™ *ultra*™ charged aerosol detector is mass sensitive and can be added to the traditional HPLC-UV or LC-MS platform. This detector provides the most consistent response for all nonvolatile and some semivolatiles of all HPLC detection techniques.¹ The detector works by charging particles (see Figure 1) and is not dependent on light scattering which can vary between analytes. The work presented here examines the combination of UV and MS detection already employed in this field, with charged aerosol detection for quantification of drug metabolites in two specific cases. This work was completed in collaboration with Bristol-Myers Squibb (BMS), using a mix of HPLC equipment and method conditions.² Although one analytical system and condition is emphasized, a variety of analytical equipment was used throughout the study.

FIGURE 1. Charged aerosol detector flow path schematic.



26802

Methods

Sample Preparation

The buspirone and erythromycin standards were incubated at 1, 30, 60, and 100 μM substrate concentrations of human liver microsome (HLM). At the end of the incubation period, the samples were quenched with an equal part of acetonitrile. The final analytical concentration of the metabolite samples discussed in this work is therefore one half of the concentration listed as the sample name (i.e., 60 μM pre-equals 30 μM postincubation).

Analytical Conditions

Column: SB C18, 1.8 μm , 4.6 \times 150 mm
Mobile Phases: A: 0.1% Formic acid in water,
B: 0.1% Formic acid in acetonitrile
Flow Rate: 1 mL/min
Total Run Time: 15 min

Detection Conditions

Thermo Scientific Dionex UltiMate™ 3000 Diode Array Detector (Thermo Scientific Accela PDA detector used in some examples):

UV Wavelength 1: 220 or 210 nm depending on example (values listed)
UV Wavelength 2: 254 nm

Dionex Corona *ultra*:

Filter: High
Gas: 35 psi Nitrogen
Range: 100 pA full scale

LTQ Orbitrap high-resolution mass spectrometer

Full scan MS with 1 \times DDS

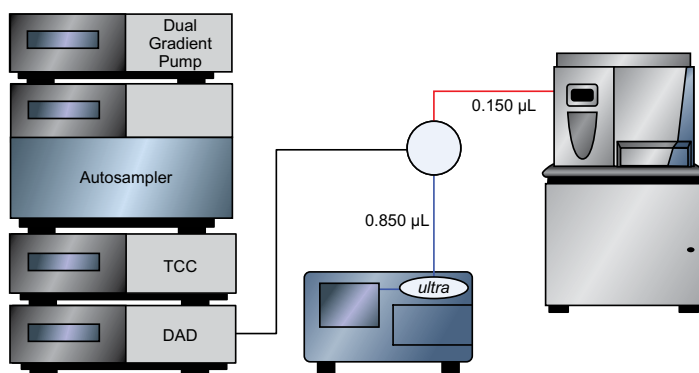
Flow Split

Valco three-way splitter

Volumes fixed by backpressure of different tubing diameters and lengths

Volume measured at charged aerosol detector inlet.

FIGURE 2. Schematic of one of the LC-charged aerosol detector-MS systems used for analysis, with a total flow of 1 mL/min split ~5.7/1 to the charged aerosol detector.



28120

Results

Buspirone

FIGURE 3. Structure for buspirone and five of its known metabolites.³

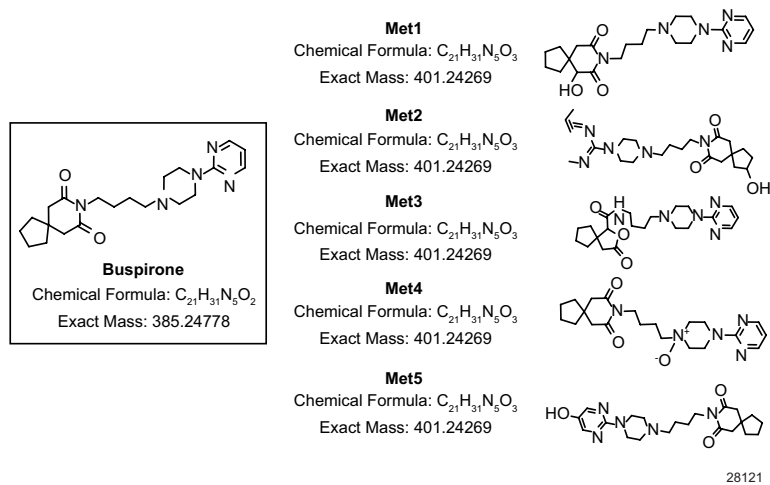


FIGURE 4. Chromatographic results for 60 μ M buspirone before and after 120 min incubation in HLM.

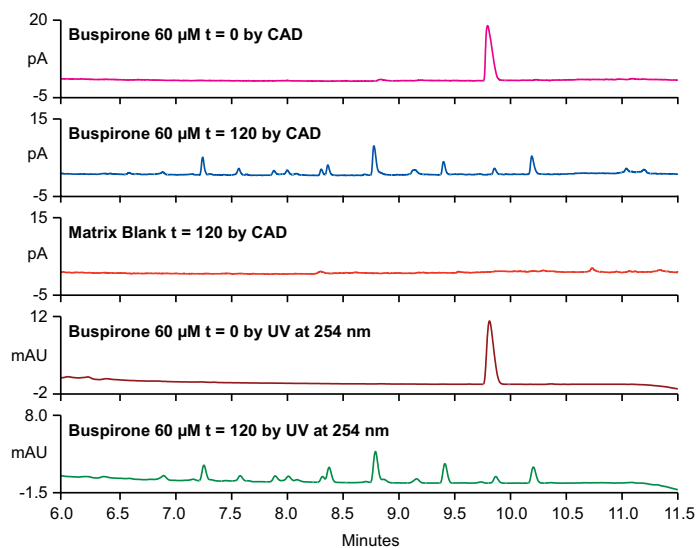


FIGURE 5. Response as peak area for injection of equal concentrations of buspirone and the five metabolites. Response deviation of each detection technique is calculated as percent RSD.

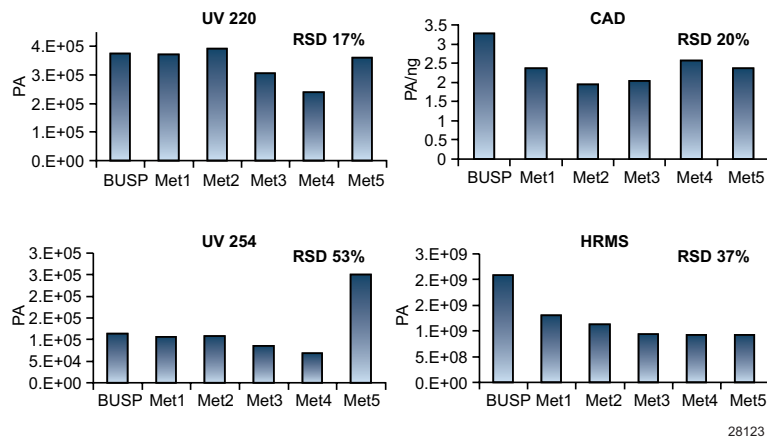
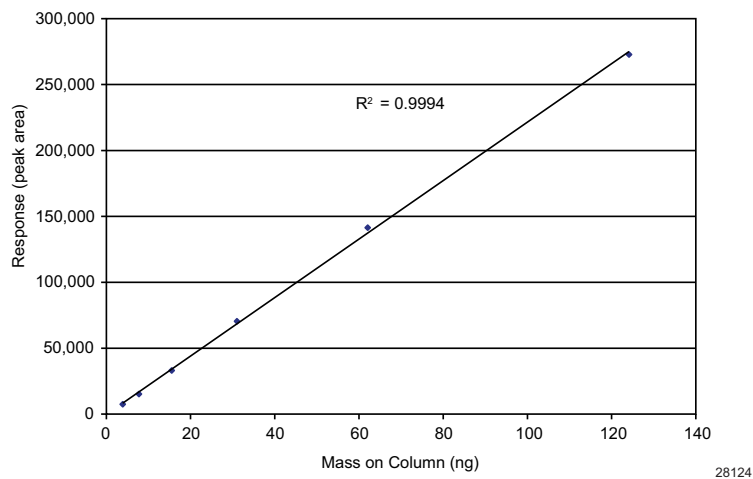


FIGURE 6. Linear response curve for buspirone with charged aerosol detection from ~4 to 124 ng on column.



Erythromycin

FIGURE 7. Results for charged aerosol detection pre- and postincubation and UV at 210 nm preincubation the 100 μM erythromycin sample.

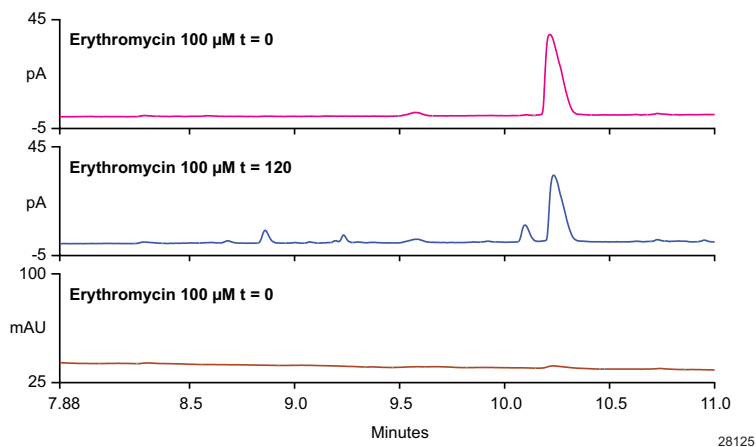
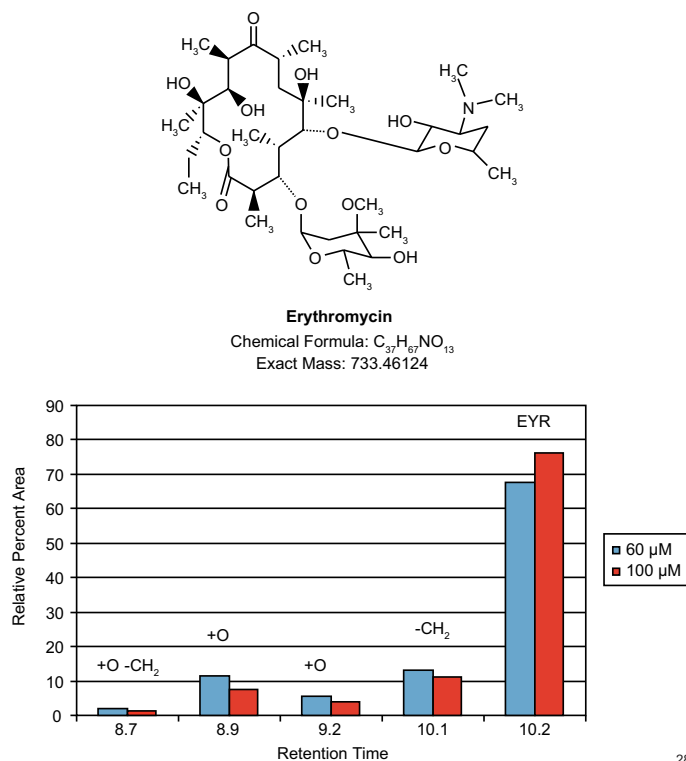


FIGURE 8. Structure (top) of erythromycin. Plot of relative area response of the four metabolites and the parent peak with charged aerosol detection (bottom).



Discussion

The results discussed here were obtained in collaboration with BMS. The optimization of the system and development of the methods were ongoing throughout the work. Several components were identified as crucial to maintaining quality data on both the charged aerosol detector and mass spectrometer. The column diameter and load volume was important, as peak splitting and shifting was observed with injection volumes greater than 30 μ L. The configuration of the flow splitting when running the LC-charged aerosol detector-MS system was determined to be the most crucial factor. The flow splitter must be positioned close to the mass spectrometer inlet using 0.004 inch or smaller i.d. tubing from the splitter to the mass spectrometer (Figure 2). The optimal flow-split range is estimated between 100 to 200 μ L/min flowing to the mass spectrometer, and the remaining flow to the charged aerosol detector. Because of the levels of sensitivity required for these analyses, a minimum flow rate of 500 μ L/min is recommended for the charged aerosol detector.

The first example, buspirone, was chosen because it is a well-characterized example, has a strong UV chromophore, and the standard material is readily available. The post HLM incubation of buspirone for both the charged aerosol detector and UV (Figure 4) show a similar metabolite profile. The five major metabolites shown in Figure 3 were quantified as relative peak areas from the initial 60 μ M buspirone sample. A sixth, minor metabolite was a volatile fragment of the buspirone and was not detected using the charged aerosol detector. The standards for the buspirone and the five metabolites were prepared and analyzed at concentrations of 15 μ M. The results for the two UV wavelengths (charged aerosol detection and MS) are shown in Figure 5. The greatest variability was found in the UV trace at 254 nm with an RSD of 53% followed by the MS at 37%. The charged aerosol detector and low wavelength UV had similar deviations of 20 and 17%, respectively.

The charged aerosol detector is typically linear from its limit of quantification to ~500 ng on column, although not a linear detector over the full dynamic range. In Figure 6, the buspirone standard was analyzed using a UHPLC method from 3.9 to 124 ng on column. The six-point calibration curve exhibited excellent linearity over this range with a correlation coefficient of 0.9994. The concentration range of interest needed for the testing of drug metabolites typically falls in this mass-on-column range, which is linear for the charged aerosol detector independent of analyte.

The second example, erythromycin, was chosen because it is representative of an array of drug candidates that fall into the nonvolatile analyte category with weak or no UV chromophores. As shown in Figure 7, the preincubation 100 μM standard has virtually no response with low wavelength UV. After incubation, the charged aerosol detector was able to detect the four major metabolites observed by MS. The relative area response for those four metabolites and the parent peak were evaluated at the 100 and 60 μM levels (Figure 8). As expected, the relative intensities of the metabolites increased at the lower concentration over the same incubation time.

The level of detection on the charged aerosol detector for the buspirone was determined to be ~ 1 ng on column with the short run UHPLC chemistry. This is typical for nonvolatile analytes under optimized conditions. Depending on molecular weight of the compound and the injection volume used, the molar concentration limit of detection can vary. While more sensitive than other nebulizer-based detection techniques, charged aerosol detection still may not provide sufficient sensitivity to meet the detection levels needed for in vivo studies. The MetPro practice (advocated by J. Joseph in 2009), of normalizing the MS metabolite ratios at a single, higher-level point with UV data then transferring that to the MS results at the low levels has been employed in previous work.¹ This was examined with the charged aerosol detector results during the calibration and is further described in Cai, et al, (2010).² This offers a method to reduce the analytical deviations due to ionization variability, while keeping the sensitivity offered by the mass spectrometer.

Conclusion

The UHPLC-UV-charged aerosol detector-MS system described in this study was able to provide crucial information needed to meet requirements of the 2008 FDA MIST initiative. The use of charged aerosol detection as confirmation of low-wavelength UV quantification for chromophoric compounds can provide additional confidence to the results. In areas where more specific wavelengths are used, such as the buspirone 254 nm, the charged aerosol detector can highlight large under- or overestimation as shown with metabolite.³ The area where this technology is most applicable is for parent or drug compounds that do not contain a strong UV chromophore. In these cases, laboratories are often left with no other option but to accept the MS quantification during the early drug discovery stages. This can lead to major issues in later drug development stages. The charged aerosol detector—while not a stand-alone solution— can provide those additional pieces of data to make accurate interpretations of in vitro data without excessive cost or time requirements.

Acknowledgments

The authors would like to thank Hong Cai and colleagues at Bristol-Myers Squibb for their work on this collaboration. Generation of the data used would not have been possible without their help.

References

1. Dependence of Response on Chemical Structure, ESA—A Dionex Company, Application Brief 70-8913 Rev B.
2. Cai, et. al. *A Novel Detection Technology: Charged Aerosol Detection (CAD) Coupled with HPLC, UV, and LTQ-Orbitrap MS for Semi-Quantitation of Metabolites in Drug Discovery Metabolism Studies*, ASMS oral presentation, Salt Lake City, UT, May, 2010.
3. Zhu. et al. *Drug Metab. Dispos.* **2005**, 33, 500–507.

www.thermoscientific.com/dionex



Thermo Scientific Dionex products are designed, developed, and manufactured under an ISO 9001 Quality System.

©2011 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

U.S./Canada	(847) 295 7500	Denmark	(45) 36 36 90 90	Sweden	(46) 8 473 3380	India	(91) 22 2764 2735
Brazil	(55) 11 3731 5140	France	(33) 1 39 30 01 10	Switzerland	(41) 62 205 9966	Japan	(81) 6 6885 1213
Austria	(43) 1 616 51 25	Germany	(49) 6126 991 0	United Kingdom	(44) 1276 691722	Korea	(82) 2 2653 2580
Benelux	(31) 20 683 9768	Ireland	(353) 1 644 0064	Australia	(61) 2 9420 5233	Singapore	(65) 6289 1190
	(32) 3 353 42 94	Italy	(39) 02 51 62 1267	China	(852) 2428 3282	Taiwan	(886) 2 8751 6655

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific