

# Simultaneous Determination of Metformin and its Chloride Counterion Using Multi-Mode Liquid Chromatography with Charged Aerosol Detection

Xiaodong Liu and Mark Tracy, Thermo Fisher Scientific, Sunnyvale, CA, USA

## Key Words

API, Counterion, Acclaim Trinity P2, Corona Veo charged aerosol detector

## Abstract

This work demonstrates the determination of metformin (API) and its counterion (chloride) using a simple isocratic method. Thermo Scientific™ Acclaim™ Trinity™ P2 columns offer cation-exchange, anion-exchange and HILIC retention mechanisms on the same phase, thus they are ideal for separating both cationic and anionic species simultaneously.

## Introduction

Determinations of Active Pharmaceutical Ingredients (API) and counterions are important assays in pharmaceutical analysis. Due to the difference in charge and/or hydrophobicity, APIs and counterions are usually analyzed by different chromatographic methods that require different separation columns and/or different instrumentation platforms. Metformin is an oral anti-diabetic drug, often formulated in its hydrogen chloride salt form. Because of the highly hydrophilic nature of both API and counterion, it is impossible to assay both components within the same analysis on any RP, ion-exchange or HILIC column.

The Acclaim Trinity P2 column is based on Nanopolymer Silica Hybrid (NSH™) technology. It consists of high-purity porous spherical silica particles coated with charged nanopolymer particles: the inner-pore area of the silica particles is modified with a covalently bonded hydrophilic layer that provides cation exchange retention while the outer surface is modified with anion-exchange nano-polymer beads. This chemistry design ensures spatial separation of the anion exchange and cation exchange regions.



Due to the fact that chloride cannot be detected by UV and that both metformin and chloride are non-volatile, an aerosol based detector is ideal for this application. The Thermo Scientific™ Dionex™ Corona™ Veo Charged Aerosol Detector represents the latest advancement of this technology. Compared with other aerosol based detectors (e.g., ELSD), the Corona Veo features superior limit of detection, better reproducibility (RSD), and ease of use. When combined with UV detection, Corona Veo is the detector of choice for many applications, including simultaneous separation of API and counterion.

The work here describes methods for determination of metformin (API) and chloride (counterion) using a simple isocratic method.

## Experimental Details

Consumables	Part Number
Acetonitrile, Fisher Optima™ LC/MS grade	A955
Formic acid, >98%	
Ammonium formate, 99.995%	
Deionized water	
Metformin hydrochloride	

Separation Conditions	Part Number
Instrumentation:	Thermo Scientific™ Dionex™ UltiMate™ 3000 RS system
Column:	Acclaim Trinity P2, 3 μm, 50 × 3 mm 085433
Mobile phase A:	Acetonitrile
Mobile phase B:	100 mM ammonium formate, pH 3.65 (6.35 g/L NH <sub>4</sub> HCO <sub>2</sub> + 4.5 g/L HCO <sub>2</sub> H)
Isocratic elution:	80% A, 20% B (v/v)
Flow rate:	0.50 mL/min
Column temperature:	30 °C
Injection volume:	1 μL
Detection:	Corona Veo Charged Aerosol Detector (evaporator temperature 55 °C, gas pressure 60 psi, data rate 5 Hz, filter 2 s, power function 1.50)

## Data processing

Software:	Thermo Scientific™ Dionex™ Chromeleon™ 6.8 SR13
-----------	---

## Results

Since Acclaim Trinity P2 columns provide both cation-exchange and anion-exchange retention mechanisms at the same time, they can adequately retain both metformin (cationic) and chloride (anionic) under the same chromatographic conditions. The unique chemistry of the Acclaim Trinity P2 column, in which cation-exchange and anion-exchange regions are spatially separated, allows for great flexibility in method optimization by adjusting mobile phase buffer concentration, pH, and/or organic solvent content. To optimize this particular application, various buffer concentrations and solvent levels were examined by proportioning acetonitrile, 100 mM ammonium formate buffer and de-ionized water. Several separation conditions were developed. The best result, according to the criteria of retention ( $k > 2$ ), resolution ( $R_s > 2$ ) and analysis time (<10 min), was achieved at 80% acetonitrile and 20% buffer, demonstrated in Figure 1.

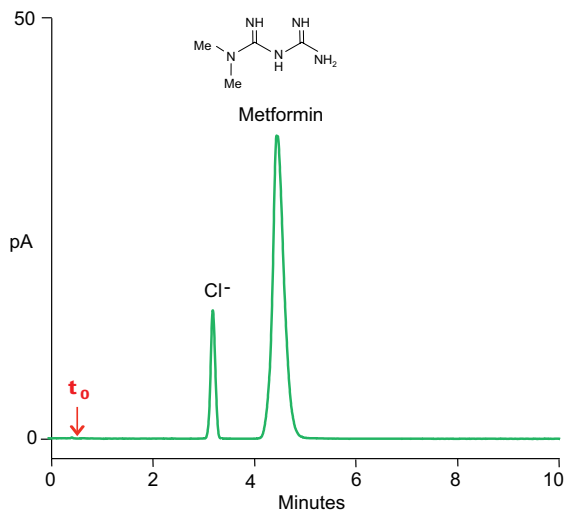


Figure 1: Simultaneous Separation of Metformin (API) and Chloride (Counterion)

## Conclusion

- The Acclaim Trinity P2 column provides solutions for simultaneous separation of API (metformin) and counterion (chloride)
- The separation is carried out using a simple mobile phase system of acetonitrile and ammonium formate buffer

## References

- [1] P.H. Stahl, C.G. Wermuth, Handbook of Pharmaceutical Salts; Properties, Selection, and Use, VCH and Wiley-VCH, Zurich, Switzerland, and Weinheim, Germany, 2008.
- [2] U.S. Food and Drug Administration Orange Book, <http://www.fda.gov/cder/ob/>.

[thermoscientific.com/specialtyLCcolumns](http://thermoscientific.com/specialtyLCcolumns)

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

**USA and Canada** +1 800 332 3331  
**Australia** 1300 735 292 (free call domestic)  
**China** 800 810 5118 (free call domestic)  
400 650 5118  
**France** +33 (0)1 60 92 48 34  
**Germany** +49 (0) 2423 9431 20 or 21

**India** +91 22 6742 9494  
+91 27 1766 2352  
**Japan** 0120 753 670 (free call domestic)  
0120 753 671 fax  
**United Kingdom** +44 (0) 1928 534 110  
**New Zealand** 0800 933 966 (free call domestic)  
**Singapore** +65 6289 1190  
**All Other Enquiries** +44 (0) 1928 534 050

**Technical Support**  
**For advice and support,**  
**please visit our website:**  
[www.thermoscientific.com/chromexpert](http://www.thermoscientific.com/chromexpert)

**Thermo**  
SCIENTIFIC

Part of Thermo Fisher Scientific