# **LC-MS Analysis of Reserpine Using High pH Buffer**

### ABSTRACT

Reserpine is being used as one of the gold standards for calibration and sensitivity testing in liquid chromatography coupled to mass spectrometry. In most cases, chromatography is performed using methanol and acetic acid (or formic acid) to improve the ionization of reserpine. A high pH buffer can be used for the LC-MS analysis of reserpine. Detection of trace level reserpine by LC-MS is now possible using high pH mobile phase.

### **OVERVIEW**

- Purpose: analysis of reserpine using a high pH buffer and comparison of the peak response (area, height and peak shape) with conventional acid containing mobile phase was performed.
- Methods: peak responses were monitored by selected ion monitoring (SIM) using a single quadrupole mass spectrometer; a parallel reaction monitoring method was established using high resolution Q Exactive<sup>™</sup> Hybrid Quadrupole-Orbitrap<sup>™</sup> mass spectrometer.
- Results: superior peak shape and higher peak intensity for reserpine was obtained with high pH buffer under the analysis conditions

### INTRODUCTION

- Reserpine is used as a calibration standard for evaluating the performance of mass spectrometers (MS) such as Thermo Scientific ion trap and triple quadrupole MS. Addition of formic acid or acetic acid to the mobile phase is necessary for sufficient ionization of reserpine. We report the use of a high pH buffer to monitor the signal response of reserpine by SIM in single quadrupole MS and compared it to the response obtained using acidic mobile phase.
- UHPLC system coupled with MS enables the generation of narrow analyte peaks in chromatographic elution. High resolution MS operated in SRM or multiple reaction monitoring (MRM) enables an acquisition frequency that is well suited for these narrow peaks generated by UHPLC.
- Reserpine was analyzed with high resolution mass spectrometry. Peak response was monitored using three different columns; both acidic and basic mobile phases were used to compare the peak response by parallel reaction monitoring. A superior peak response was obtained using high pH buffer.

## METHODS

Instrumentation

- Thermo Scientific high resolution Q Exactive Hybrid Quadrupole-Orbitrap mass spectrometer with Vanquish™ UHPLC system
- Agilent 1260 system and 6150 MS (single quadrupole) with jet stream ion focusing technology, diode array detector (DAD) and auto-sampler
- ESI-MS: Agilent 1100 with MSD SL (single quadrupole)
- Columns:
- Agilent Poroshell 120 EC-C18 column (50 x 4.6 mm, 2.7 micron) - Agilent XDB extend C18 column (50 x 2.1 mm, 3.5 micron)
- Thermo Scientific Acclaim<sup>™</sup> 120 column (50 x 3.0 mm, 3 micron)
- Mobile phases and reagents:

- Water (W8), methanol (A458), formic acid (A117), acetonitrile (A956), ammonium bicarbonate, and ammonium hydroxide solution are obtained from Fisher Scientific (Pittsburgh, PA)

- Reserpine was purchased from Sigma Chemical Company (St. Louis, MO) • LC-MS testing was performed using three different mobile phases:
- 10 mM ammonium bicarbonate, pH 10 with 5% methanol as mobile phase A and methanol as mobile phase B
- 0.1% formic acid in water and acetonitrile - 0.1% acetic acid in water and methanol
- LC-MS run conditions:
- Flow rate 0.5 mL/min and 0.3 mL/min; injection volume 5  $\mu$ L
- Gradient runs were performed as follows:

### I C with single quadrupole MS

Lo mai single quadrupole mo					
Time (min)	Mobile phase A%	Mobile phase B%	Time (min)	Mobile phase A%	Mobile phase B%
0	60	40	0	60	40
2	60	40	2	60	40
10	0	100	5	0	100
20	0	100	7	0	100

• Mass spectrometer (single quadrupole) data acquisition condition:

- Fragmentor: 70 V
- Dry gas temperature: 250° C
- Dry gas flow: 12 mL/min
- Capillary voltage: 3000 √ - Gain: 1
- Threshold: 150
- Dwell msec: 290
- Peak width: 0.1 min
- Cycle time: 0.84 sec/cycle
- High resolution hybrid mass spectrometer data acquisition condition:
- Full MS SIM
- Polarity: Positive - Resolution: 70,000
- Microscans: 1
- AGC target: 1e6
- Scan ranges: 50 750 m/z
- Maximum IT: 200 msec
- CE: 40

LC	LC with high resolution MS		
	Time (min)	Mobile phase A%	Mobile phase B%
	0	60	40
	2	60	40
	5	0	100

with	high	resolution	MS



-igure 2. Linear Res		
PPB Conc.	Are	
1	1.	
2.5	2.	
10	7.	
25	1.	
50	3.	
100	5.	



Conc (ppb)	Res_Ar
1	6.0
2.5	1.1
25	1.2
50	3.5
100	4.1
200	7.0

Subhra Bhattacharya and Stephen C. Roemer, Thermo Fisher Scientific, Laboratory Chemicals, One Reagent Lane, Fair Lawn, NJ 07410