

# Extraction of Basic Drugs from Plasma with Polymeric SPE

## Application Note

Pharmaceuticals

### Authors

William Hudson, David Jones,  
Arnie Aistars, and Max Erwine  
Agilent Technologies, Inc.  
5301 Stevens Creek Boulevard  
Santa Clara, CA 95051  
USA

### Introduction

Bioanalytical solid phase extraction (SPE) has been dominated by polymeric sorbents in recent years. The ease-of-use, good flow, and resistance to effects of drying relative to silica-based sorbents make polymeric sorbents an obvious choice for high volume, high throughput assays requiring quick validation and minimal method development.

Because the method validation process is time consuming and requires high quality data, SPE methods that are fast, yet produce good recoveries with high reproducibility, are desirable. To the extent that the SPE process is streamlined without compromising data integrity, method validation can be simplified and shortened. Bond Elut Plexa minimizes method development with simple and effective methods and improves analytical sensitivity and reproducibility with an advanced polymeric structure that minimizes binding of large biomolecules to the surface, with the end result of simplifying and streamlining the SPE process.



**Agilent Technologies**

## Materials and Methods

### SPE reagents and solutions

2% ammonium hydroxide    Add 20  $\mu$ L concentrated ammonium hydroxide to 1 mL DI H<sub>2</sub>O

Methanol                Reagent grade or better

5% methanol            Add 5 mL methanol to 95 mL DI H<sub>2</sub>O

Bond Elut Plexa        10 mg 96 well plate (p/n A4969010)

### SPE method

Sample                 100  $\mu$ L human plasma

Pretreat                Dilute with 300  $\mu$ L 2% NH<sub>4</sub>OH

Condition              1. 500  $\mu$ L CH<sub>3</sub>OH  
                              2. 500  $\mu$ L H<sub>2</sub>O

Wash                    500  $\mu$ L 5% CH<sub>3</sub>OH in H<sub>2</sub>O

Elute                    500  $\mu$ L CH<sub>3</sub>OH

All samples evaporated to dryness and reconstituted in 100  $\mu$ L of 80:20 0.1% formic acid: CH<sub>3</sub>OH aq.

LC/MS performed – ESI, drying gas @ 400 °C, 30 psi

## LC conditions

### Mobile phase

A                        0.1% Formic acid

B                        Methanol

### LC gradient program

Time (min)    %B

0:00                40

0:15                40

1:00                80

3:00                80

4:30                40

### Column

Type                 Pursuit XRs C18 3  $\mu$ m, 50  $\times$  2.0 mm (p/n A3001050X020)

Flow rate            0.2 mL/min

## Results and Discussion

The procedure described provides a simple and effective SPE method for the extraction of basic or neutral drugs from human plasma. The Limit of Quantitation (LOQ) of the combined SPE and LC/MS/MS analysis was 1.0 ng/mL. The internal standard for the application was 50 ng/mL quetiapine.

Recoveries were calculated from a second order regression with RSD values based on a sampling of n = 6. Excellent recoveries were achieved demonstrating good retention and elution, as well as minimal ion suppression. Response for all the compounds evaluated was linear up to three orders of magnitude from 1.0 ng/mL to 1.0  $\mu$ g/mL with correlation coefficients all above 0.995 (n = 6). To demonstrate reproducibility, samples were analyzed at two concentrations (n = 6). Figure 1 shows the chromatograms of the extractions at 100 ng/mL. As shown in Table 1, the extractions produced reproducibly high recoveries.

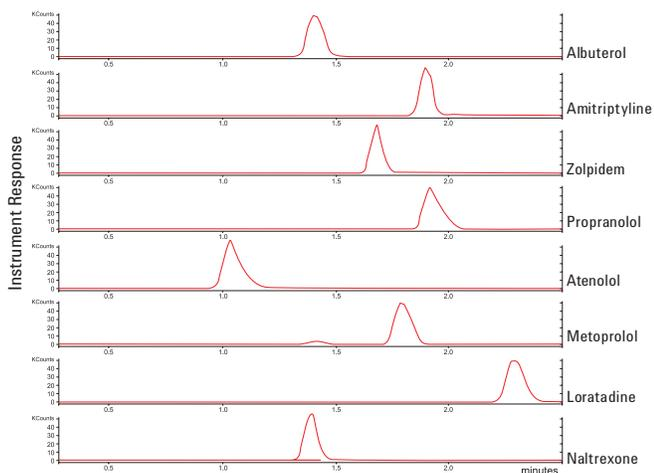


Figure 1. Chromatograms of a 100 ng/mL human plasma extract.

Table 1. High Recoveries of Basic Drugs with Bond Elut Plexa

Drug	log P	pKa	0.5 µg/mL %Recovery	%RSD	1.0 µg/mL %Recovery	%RSD
Albuterol	1.3	10.3	95	5	100	2
Amitriptyline	4.6	9.4	100	10	100	4
Zolpidem	3.9	6.2	100	8	103	2
Propranolol	3.6	9.5	102	6	101	6
Atenolol	1.3	9.6	97	4	101	4
Metoprolol	1.3	10.8	100	5	100	5
Loratadine	5.2	4.9	97	5	95	3
Naltrexone	1.8	9.2	103	11	100	4

## Conclusions

Bond Elut Plexa is a useful tool for high-throughput SPE applications that require analysis at low analyte levels, need validated reproducibility, and must be quickly implemented with minimal method development. A single method for basic analytes covers a broad range of analyte polarities and delivers reproducibly high recoveries. Bond Elut Plexa is therefore highly recommended for bioanalytical work in pharmaceutical clinical research trials, including contract research.

## For More Information

These data represent typical results. For more information on our products and services, visit our Web site at [www.agilent.com](http://www.agilent.com).

[www.agilent.com](http://www.agilent.com)

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

Information is subject to change without notice.

© Agilent Technologies, Inc.  
Printed in the USA  
March 21, 2011  
5990-7685EN



**Agilent Technologies**