# Customer Application Note

## Determination of the Punicalagins Found in Pomegranate by High Performance Liquid Chromatography

can 106

Steve Baugh, Jessica Revell, and Kim Eastman Chromadex, Irvine, CA 92618

## Introduction

Antioxidants found in fruits and vegetables or the juices of fruits and vegetables are known to provide our bodies with many health benefits. Pomegranate is one such fruit that contains high levels of antioxidants; studies have shown that pomegranate contains more antioxidants than green tea, cranberries, and even red wine. The antioxidants in pomegranates include polyphenols, such as tannins and anthocyanins. The most abundant polyphenols in pomegranate juice are the hydrolyzable tannins called punicalagins that have freeradical scavenging properties in laboratory experiments.

For more details on pomegranate antioxidant properties click the following link:

http://pomegranateinformation.com/pomegranate/antioxidant-properties/

## Summary

Samples			
-	Lot #	CDXA #	
Punicalagins A & B Chromadex	16983-1218	CDXA-119.113	

## Results

See chromatograms in Data section:

## Analytical Method Laboratory Supplies

Analytical Balance Ultrasonication Bath Assorted Volumetric glassware Syringes and Syringe Filters HPLC/GC glass vials and caps

## Solvents and Reagents

Formic Acid Milli-Q<sup>®</sup> Water Acetonitrile (CH<sub>2</sub>CN)



## Solution Preparation

Mobile Phase: 1 % Formic Acid in Milli-Q water The solution was prepared by adding 10 mL of Formic Acid to 1 L of Milli-Q water and mixing well.

## Stock Standard Preparation

The stock standard was prepared by weighing 5.42 mg of punicalagins A & B into a 10 mL volumetric flask. 5 mL of Milli-Q water were added and sonicated. The solution was then allowed to re-equilibrate to ambient temperature, brought to volume with Milli-Q water, and mixed well.

## Instrument Parameters

Instrument:	Dionex HPLC System		
Detection:	UV-vis		
Mobile Phase A:	1% Formic Acid in Milli-Q Water		
Mobile Phase B:	Acetonitrile		
Gradient Program:			
Time (min)	%A	%В	
0	95	5	
18	85	15	
20	35	65	
25	95	5	
30	95	5	
Column:	Dionex Acclaim® Polar Advantage PA II		
	150 × 3.0	0 mm, 3 µm particle	
Flow Rate:	1.0 mL/min		
Temperature:	30 °C		
Injection Volume:	10 µL		
UV Detection:	260 nm		



## Data

**Figures** 

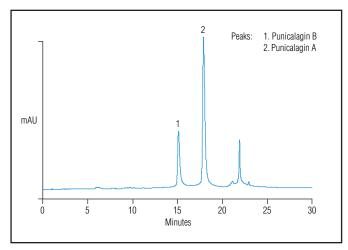


Figure 1. Punicalagins mixed standard injection.

## Structure

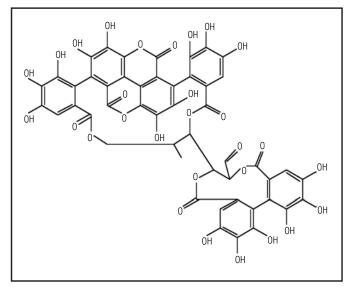


Figure 2. Punicalagin.

## References

- 1. ChromaDex Analytics Laboratory Notebook 73, pages 114–115.
- 2. ChromaDex Analytics Laboratory Notebook 119, page 113.
- 3. ChromaDex SOP "Routine Laboratory Calculations."
- 4. Analytical Method: 99.1-CD-1.0-000216.

Milli-Q is a registered trademark of Millipore. Acclaim is a registered trademark of Dionex Corporation.

## Passion. Power. Productivity.

## DIONEX

## **Dionex Corporation**

1228 Titan Way P.O. Box 3603 Sunnyvale, CA 94088-3603 (408) 737-0700

#### North America

Europe

#### U.S./Canada (847) 295-7500

**South America** Brazil (55) 11 3731 5140 Austria (43) 1 616 51 25 Benelux (31) 20 683 9768; (32) 3 353 4294 Denmark (45) 36 36 90 90 France (33) 1 39 30 01 10 Germany (49) 6126 991 0 Ireland (353) 1 644 0064 Italy (39) 02 51 62 1267 Sweden (46) 8 473 3380 Switzerland (41) 62 205 9966 United Kingdom (44) 1276 691722

#### Asia Pacific

Australia (61) 2 9420 5233 China (852) 2428 3282 India (91) 22 2764 2735 Japan (81) 6 6885 1213 Korea (82) 2 2653 2580 Singapore (65) 6289 1190 Taiwan (886) 2 8751 6655

**ISO** 9001

www.dionex.com

