

Fast Analysis of Triterpene Glycosides in Black Cohosh Using the Agilent 1290 Infinity LC System and an Agilent Poroshell 120 SB C-18 2.7 μm Column

Application Note

Author

Nick Severin
Schwabe North America,
Green Bay, WI
USA

Abstract

Using the Agilent 1290 Infinity LC System coupled with the Agilent Poroshell 120 SB C-18, 3 \times 100 mm, 2.7 μm column can not only reduce solvent consumption, but allows for greater peak resolution as well. This is illustrated by comparing analyses of triterpene glycosides in black cohosh extract using the parameters outlined in the United States Pharmacopeia (USP) and the 1290 Infinity LC System with a Poroshell 120 SB C-18 column.



Agilent Technologies

Black Cohosh Supplements

Black Cohosh (*Actaea racemosa* and *Cimicifuga racemosa*) is a member of the buttercup family, native to North America. Extracts of black cohosh are made from its roots and rhizomes. The extracts are standardized to the amount of triterpene glycoside content using the 26-deoxyactein standard. Black cohosh extract is mainly used to treat menopausal symptoms. It is regulated as a dietary supplement, rather than as a drug.

Black Cohosh Extraction Conditions

Black cohosh extract source Martin Bauer Group, Germany

Lot used 11018338

A 300 mg amount of sample extract was dissolved in 10 mL of methanol with the aid of a 20 minute sonication step.

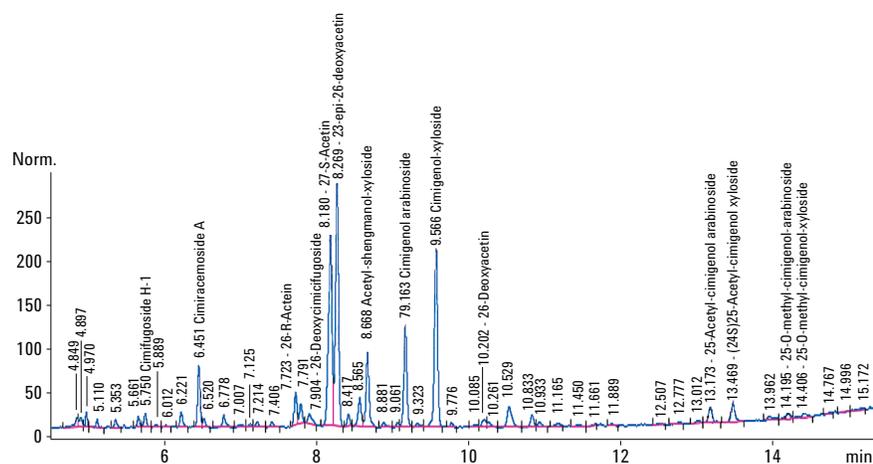


Figure 1A. Triterpene glycosides in black cohosh extract using an Agilent 1290 Infinity LC System with an Agilent Poroshell 120 SB C-18, 2.7 μ m column.

Instrument parameters for Figures 1A and 1B

Column Agilent Poroshell 12,0 SB C18, 3 \times 100 mm, 2.7 μ m, 35 $^{\circ}$ C

Flow rate 1.0 mL/minute

Injection volume 5.0 μ L

ELSD detection conditions

Instrument Agilent 385 Evaporative Light Scattering Detector, Model G4261A

Nebulizer temperature 40 $^{\circ}$ C

Evaporator temperature 40 $^{\circ}$ C

Nitrogen gas flow rate 1.50 SLM

PMT gain 2.0

Data rate 40 Hz

Mobile phase A 0.1% Formic acid in water

Mobile phase B Acetonitrile

Gradient

Time (min)	% A	% B
0	80	20
2.19	80	20
4.10	68	32
15.03	36	64
17.77	5	95
18.77	5	95
19.13	80	20
22.20	80	20

Results and Discussion

The chromatogram in Figure 1A shows a time reduction 3.83 times better than the USP conditions shown in Figure 2A, and a solvent savings of 6.1 times. The USP assay uses a 250 \times 4.6 mm, 5.0 μ m C-18 column. This is compared to a Poroshell C-18, 100 \times 3 mm, 2.7 μ m column using the 1290 Infinity LC System. Both methods use the same mobile phase, detector settings (Agilent 385-Evaporative Light Scattering Detector (ELSD)) and column temperature.

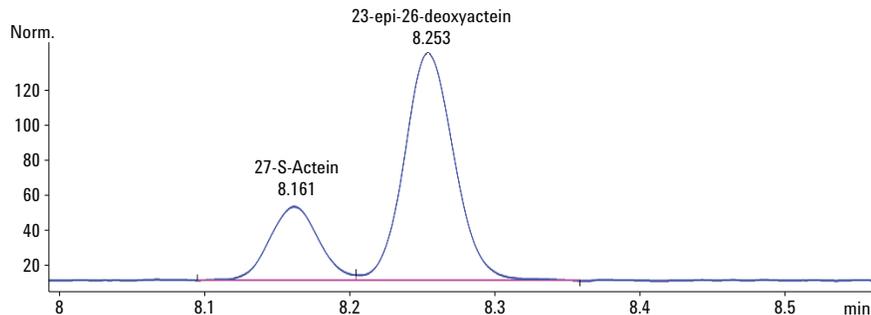


Figure 1B. Chromatogram of system suitability solution using an Agilent 1290 Infinity LC System and an Agilent Poroshell 120 SB C-18, 2.7 μ m column. Sample description: 100 μ g/mL solution of (27S) actein and 23-epi-26-deoxyactein standards.

Resolution factor 1.47
Acceptance criteria NLT 1.0

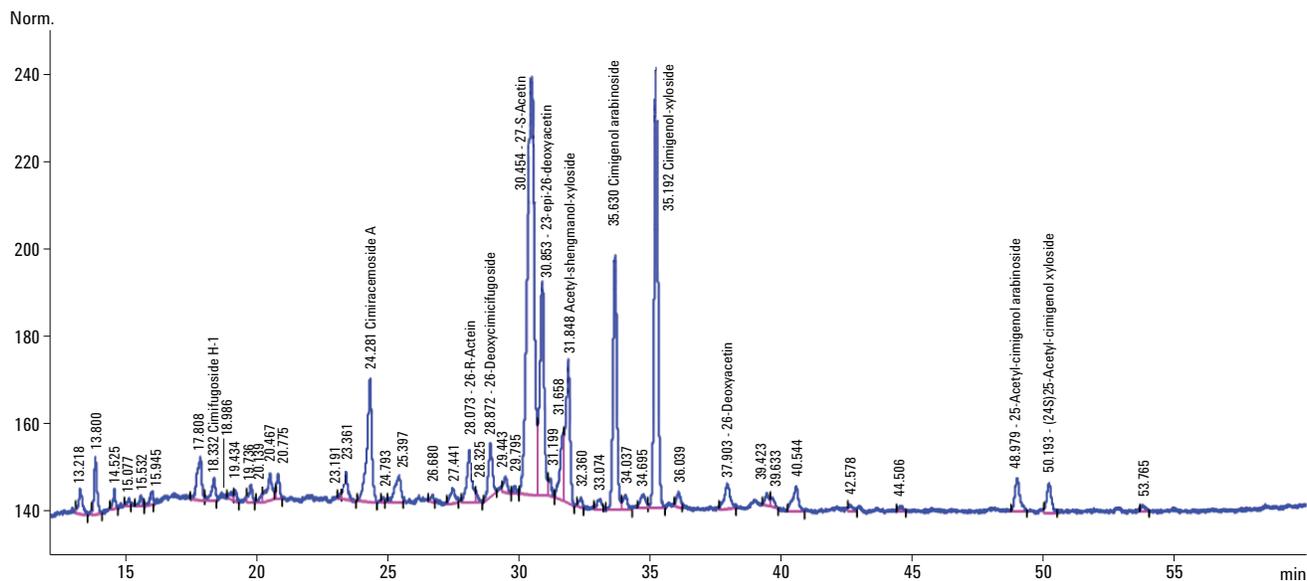


Figure 2A. Triterpene glycosides in black cohosh extract. USP method using a Phenomenex Luna C-18, 250 \times 4.6 mm, 5.0 μ m column.

Instrument parameters for Figures 2A and 2B

Column Phenomenex Luna C-18, 250 \times 4.6 mm, 5 μ m, 35 $^{\circ}$ C

Flow rate 1.6 mL/minute

Injection volume 20 μ L

ELSD detection conditions

Instrument Agilent 385 Evaporative Light Scattering Detector, Model G4261A

Nebulizer temperature 40 $^{\circ}$ C

Evaporator temperature 40 $^{\circ}$ C

Nitrogen gas flow rate 1.50 SLM

PMT gain 2.0

Data rate 40 Hz

Mobile phase A 0.1% Formic acid in water

Mobile phase B Acetonitrile

Gradient

Time (min)	% A	% B
0	80	20
8	80	20
15	68	32
55	36	64
65	5	95
70	5	95
75	80	20
85	80	20

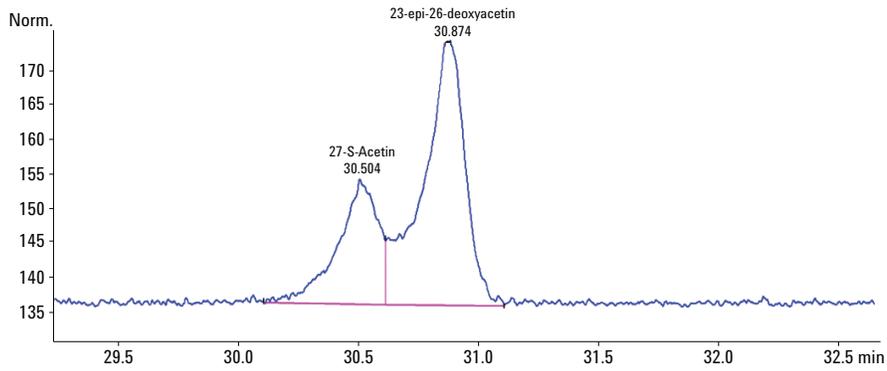


Figure 2B. Chromatogram of system suitability and a Luna 4.6 × 250 mm C18, 5µm, column. 100 µg/mL solution of (27S) acetin and 23-epi-26-deoxyacetin standards.
 Resolution factor 1.20
 Acceptance criteria NLT 1.0 USP Method

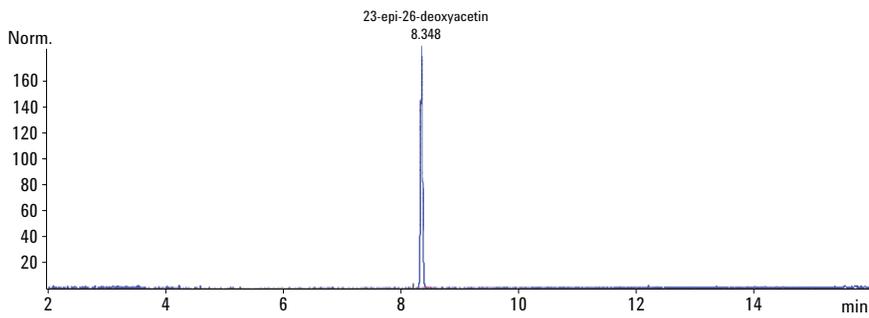


Figure 3. 23-epi-26-deoxyacetin reference standard, chromadex #00004130, Lot # 4130-512
 Calibration range 12.2–391.8 µg/mL
 R² 0.99999

Conclusion

In analyzing complex botanical extracts, the Agilent 1290 Infinity LC System, in conjunction with 3.0 mm id Agilent Poroshell 120 SB C-18, columns, can speed up analysis time, generating labor costs and solvent consumption. This is accomplished concurrently with improved resolution. The resolution factor between (27S)Actein and 23-epi-26-deoxyactein standard peaks using the USP method was 1.2 (Figure 2B) and the resolution factor for the Poroshell method was 1.47 (Figure 1B). With the current high cost of acetonitrile, a 6.1 times reduction in solvent consumption allows the analyst continue with acetonitrile in the method, avoiding the time-consuming redevelopment process, and avoid resolution changes and peak broadening noted with methanol mobile phases.

Reference

1. Black Cohosh Powder Extract Monograph, United States Pharmacopeia, USP35, NF30 p. 1210.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2013
Printed in the USA
May 2, 2013
5991-2366EN



Agilent Technologies