

Determination of Yohimbine and Related Alkaloids Using HPLC-ECD

Paul Gamache and Ian Acworth
Thermo Fisher Scientific, Chelmsford, MA, USA

Key Words

Alkaloids, HPLC-ECD, Yohimbine, Vinca

Goal

To develop a sensitive and selective HPLC-ECD method capable of measuring the low levels of yohimbine and related alkaloids typically found in plant extracts.

Introduction

Yohimbine is an indolealkylamine alkaloid with chemical similarity to reserpine. Yohimbine has been isolated from *Cornanthe johimbe*, *Rubiaceae* and related trees, as well as *Rauwolfia serpentina*. Both yohimbine and a diastereoisomer, corynanthine, are α 2-adrenoceptor antagonists of limited duration.¹ Yohimbine has limited effect on smooth muscle, but it readily passes through the blood-brain barrier and produces a complex pattern of responses at doses lower than required to produce a-adrenergic blockade. Yohimbine has been reported to produce central excitation (e.g., elevated blood pressure, increased heart rate, increased motor activity, tremor, irritability, vomiting, nausea and sweating) and to act as an anti-diuretic (through its ability to release antidiuretic hormone).² Some evidence also suggests that yohimbine can act as an aphrodisiac,³ although this effect may be limited to individuals with vascular diseases.^{4,5}

A variety of techniques have been used to measure yohimbines including: GC-MS and HPLC with UV, fluorescence or amperometric detection.⁶⁻⁸ In general these approaches suffer from a lack of sensitivity, selectivity or require extensive sample preparation and derivatization before analysis. A gradient HPLC coulometric array method was developed capable of measuring yohimbine, related alkaloids and the vinca alkaloids simultaneously.

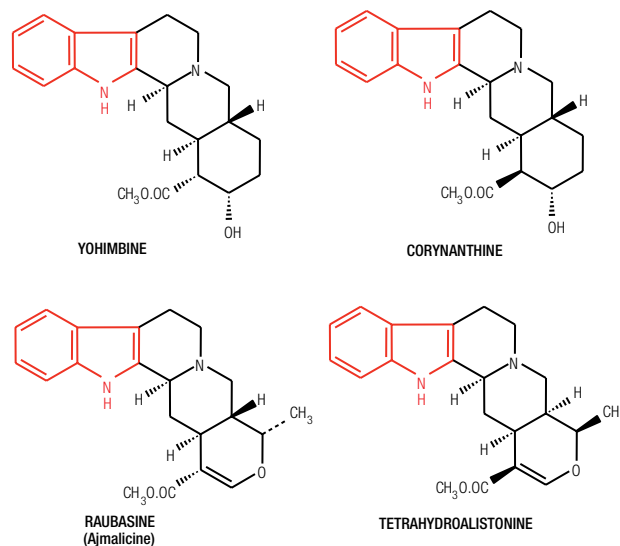


Figure 1. The structure of Yohimbine, Corynanthine, Raubasine and Tetrahydroalstonine.

This approach offers low picogram sensitivity and uses voltammetric resolution to correctly identify each analyte and check for possible coelutions. This method will have application to the measurement of these alkaloids in natural products (e.g., tree bark) and for the study of their pharmacokinetics *in vivo*.

Materials and Methods

The gradient analytical system consisted of two pumps, an autosampler, and a twelve-channel Thermo Scientific™ Dionex™ CoulArray™ Coulometric Array Detector. Higher levels of analytes were also verified using an UV absorbance detector placed after the array.

Conditions	
Column:	C18, 4.6 × 150mm, 5µm
Mobile Phase A:	100 mM Sodium acetate; acetonitrile; methanol (85:10: 5 v/v/v), final pH 6.2 with phosphoric acid
Mobile Phase B:	100 mM Sodium acetate; acetonitrile; methanol (50:30:20 v/v/v), final pH 6.2 with phosphoric acid
Gradient Conditions:	Isocratic 25%B until 1.1 minutes. Linear increase of phase B from 25% to 100% over 19 min. Hold for 18 min. Isocratic, 25% B for 3 min
Flow Rate:	1.0 mL/min
Temperature:	Ambient
Injection Volume:	20 µL
Detector Conditions	
Detector:	Model 5600A, CoulArray
Applied Potentials:	+200 to +950 mV vs. Pd. in 75 mV
Wavelength:	274 nm

Results and Discussion

Yohimbine and related alkaloids were readily resolved both chromatographically and voltammetrically from the other alkaloids in under 41 min (Figure 2). The method was linear over at least four orders of magnitude and had limits of detection of < 50 pg for all compounds examined.

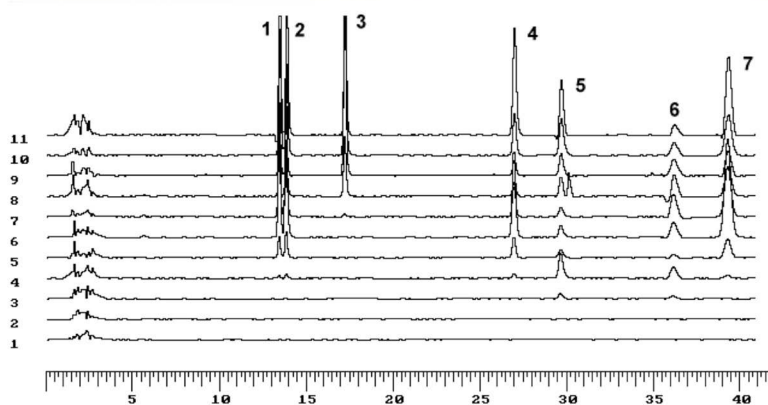


Figure 2. Chromatogram showing separation of vinca and other alkaloids (10 ng on column). 1 - corynanthine; 2 - yohimbine; 3 - vincamine; 4 - ajmalicine; 5 - vincristine; 6 - vinblastine; 7 - tetrahydro-alistinine. The applied potentials were channel 1 to channel 11, +200 to +950mV in +75mV increments. The chromatogram is presented at a gain of 500 nA.

The CoulArray detector is easily capable of measuring the low levels of yohimbine and related alkaloids typically found in plant extracts. This can be done routinely because of the inherent stability of the coulometric electrode design. As analytes are identified both by retention time and voltammetric behavior across the array, compound misidentification and coelution can be minimized.^{9,10} This approach allows analysis in complex samples such as plant extracts and biologicals with minimal sample preparation. Furthermore, this method can be used to simultaneously measure a variety of other alkaloids such as the vinca alkaloids.

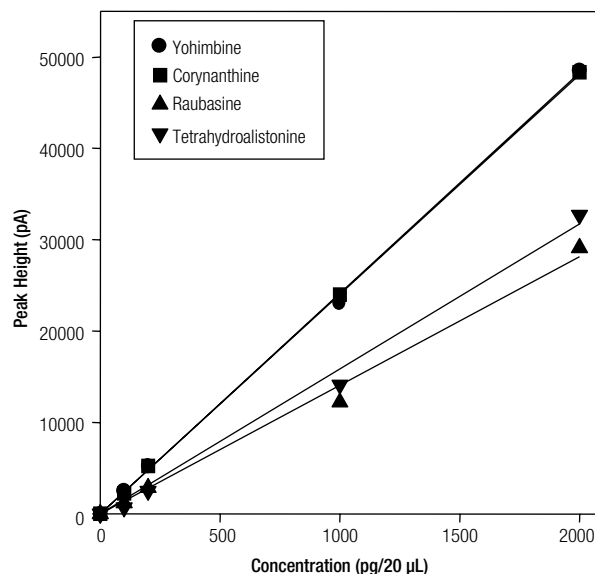


Figure 3. Linearity of yohimbine and related alkaloids.

References

- Goldberg, M. R., and Robertson, D. (1983). Yohimbine: A pharmacological probe for the alpha 2-adrenoreceptor. *Pharmacol. Rev.*, **35**, 143-180.
- Goodman Gilman, A., Goodman, L. S., and Gilman, A. (Eds.). (1980). *The Pharmacological Basis of Therapeutics*. Sixth edition. MacMillan Publishing Co., Inc. New York.
- Rosen, R. C., and Ashton, A. K. (1993). Prosexual drugs: Empirical status of "new aphrodisiacs." *Arch. Sex. Behav.*, **22**, 521-543.
- DeWire, D. M. (1996). Evaluation and treatment of erectile dysfunction. *Am. Fam. Physician*, **53**, 2101-2108.
- Montorsi, F., Guazzoni, G., Rigatti, P., and Poxxa, G. (1995). Pharmacological management of erectile dysfunction. *Drugs*, **50**, 465-79.
- Diquet, B., Doare, L., and Gaudel, G. (1984). New method for the determination of yohimbine in biological fluids by high-performance liquid chromatography with amperometric detection. *J. Chromatogr.*, **311**, 449-455.
- Reimer, G., Suarez, A., and Chui, Y. C. (1993). A liquid chromatographic procedure for the analysis of yohimbine in equine serum and urine. *J. Anal. Toxicol.*, **17**, 178-181.
- Ylinen, M., Suhonen, P., Naaranlehti, T., Lapinjoki, S., and Huhtikangas, A. (1990). Gas chromatographic-mass spectrometric analysis of major indole alkaloids of *Catharanthus roseus*. *J. Chromatogr.*, **505**, 229-234
- Acworth, I., and Gamache, P. (1996). The coulometric electrode array for use in HPLC analysis. Part 1: Theory. *American Lab.*, **5**, 33-38.
- Svendsen, C. (1993). Multi-electrode array detectors in high-performance liquid chromatography: A new dimension in electrochemical analysis. *Analyst*, **118**, 123-129

Ordering Information

Description	Part Number
HPG-3400RS Biocompatible Binary Rapid Separation Pump with two solvent selector valves	5040.0046
WPS-3000TBRS Biocompatible Rapid Separation	5841.0020
CoulArray, Model 5600A - 12 channel	70-4329
CoulArray Organizer with Temp. Control	70-4340T
Accessory Kit, CoulArray Detector to Thermo Scientific™ Dionex™ UltiMate™ 3000 System	70-9191

www.thermofisher.com/chromatography

©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 manners 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.

Africa +43 1 333 50 34 0

Australia +61 3 9757 4300

Austria +43 810 282 206

Belgium +32 53 73 42 41

Brazil +55 11 3731 5140

Canada +1 800 530 8447

China 800 810 5118 (free call domestic)
400 650 5118

Denmark +45 70 23 62 60

Europe-Other +43 1 333 50 34 0

Finland +358 9 3291 0200

France +33 1 60 92 48 00

Germany +49 6103 408 1014

India +91 22 6742 9494

Italy +39 02 950 591

Japan +81 6 6885 1213

Korea +82 2 3420 8600

Latin America +1 561 688 8700

Middle East +43 1 333 50 34 0

Netherlands +31 76 579 55 55

New Zealand +64 9 980 6700

Norway +46 8 556 468 00

Russia/CIS +43 1 333 50 34 0

Singapore +65 6289 1190

Sweden +46 8 556 468 00

Switzerland +41 61 716 77 00

Taiwan +886 2 8751 6655

UK/Ireland +44 1442 233555

USA +1 800 532 4752

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand