Determination of Vinca Alkaloids in Periwinkle Plants Using HPLC-ECD

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

Key Words

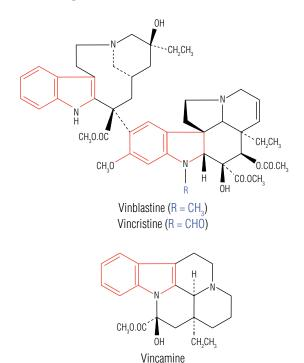
Periwinkle, HPLC-ECD, Alkaloids, Bis-Indole Derivatives, Neoplastic Diseases

Goal

To develop an HPLC-ECD method capable of measuring low levels of the vinca alkaloids typically found in plant extracts

Introduction

The vinca alkaloids consist of the bis-indole derivatives, vinblastine and vincristine isolated from the periwinkle plant, *Vinca rosea* Linn and the monomeric indole derivative vincamine, which is the major alkaloid of *Vinca minor* Linn (Figure 1). The vinca alkaloids are used as chemotherapeutic agents to treat a variety of neoplastic diseases including Hodgkin's disease, choriocarcinoma, acute and chronic leukemias, lymphosarcomas and a variety of other cancers. Their mechanism of action involves binding to tubulin with disruption of mitotic spindle formation eventually leading to inhibition of mitosis, metaphase arrest and cell death.¹





Vinca alkaloids have previously been measured using a variety of techniques including: radioassay and highperformance liquid chromatography with UV, fluoresence and electrochemical detection (ECD).^{2–5} In general these approaches suffer from either a lack of sensitivity (HPLC techniques) or selectivity (radioimunoassay and HPLC techniques). Coulometric electrode array detection when coupled to gradient HPLC possesses both the selectivity and sensitivity required to measure the vinca alkaloids in natural products (e.g. periwinkle leaf extracts) and for the study of their pharmacokinetics and metabolism in vivo. Furthermore, this method is readily adaptable for the analysis of the ever broadening variety of synthetic analogs.⁶

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.



Figure 1. The chemical structures of vinblastine, vincristine, and vincamine.

LC Conditions	
Column:	Base-deactivated, 4.6 \times 150 mm, 5 μm
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 min. 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
Detectors Conditions	
Detector:	Model 5600A, CoulArray
Applied Potentials:	+200 to +950 mV vs Pd. in 75 mV
Detector Wavelength:	274 nm

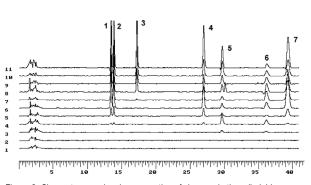
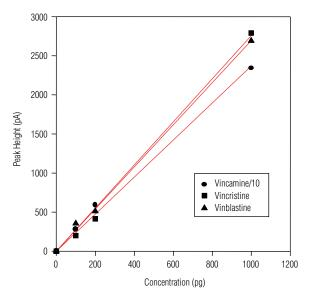
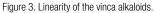


Figure 2. Chromatogram showing separation of vinca and other alkaloids (10 ng on column). 1-corynanthine; 2-yohimbine; 3-vincamine; 4-ajmalicine; 5-vincristine; 6-vinblastine; 7-tetrahydro-alistonine. The applied potentials were channel 1 to channel 11, +200 to +950 mV in +75 mV increments. The chromatogram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





Results and Discussion

The vinca 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 3).

The CoulArray detector is easily capable of measuring the low levels of the vinca 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.^{7,8} Furthermore, this method has greater sensitivity and selectivity than PDA-based approaches. This allows analysis in complex samples such as plant extracts and biologicals with a minimum of sample preparation. This same method can be used to measure a variety of other alkaloids such as yohimbine.

References

- 1. Castle, M. C. (1989). Plant alkaloids: The vinca alkaloids. Cancer Growth Prog., 10, 147–151.
- Solere, P. and Lucas, C. (1991). Analytical limits in clinical pharmacokinetics: example of vinca alkaloids. Bull. Cancer, 78, 775–787.
- Van Tellingen, O., Beijnen, J.H., and Nooyen, W. J. (1991). Analytical methods for the determination of vinca alkaloids in biological specimens: a survey of the literature. J. Pharm. Biomed. Anal., 9, 1077–1082.
- Vendrig, D. E. M. M., Teeuwsen, J., and Holthuis, J. J. M. (1988a). Analysis of vinca alkaloids in plasma and urine using high-performance liquid chromatography with electrochemical detection. J. Chromatogr. 424, 83–94
- Vendrig, D. E. M. M., Teeuwsen, J., and Holthuis, J. J. M. (1988b). Analysis of vinca alkaloids in plasma and urine using high-performance liquid chromatography with electrochemical detection. J. Chromatogr. 434, 145–155
- 6. Budman, D. R. (1992). New vinca alkaloids and related compounds. Semin. Oncol., 19, 639–645.
- Acworth, I. N., and Gamache, P. (1996). The coulometric electrode array for use in HPLC analysis. Part 1: Theory. American Lab., 5. 33–38.
- 8. Bowman, R. N. (1996). Analysis of anti-cancer alkaloids vincristine and vinblastine. Chromatography Update, Rainin.

Ordering Information

In the U.S., call (800) 346-6390 or contact the Thermo Fisher Scientific Regional Office nearest you. Outside the U.S., order through your local Thermo Fisher Scientific office or distributor. Refer to the following part numbers.

Product 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 (ree call domestic) 400 650 5118 AB71166-EN 07165 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



A Thermo Fisher Scientific Brand