



- ASE Features and Benefits
- Application Notes
- Literature References
- Customer References

### Accelerated Solvent Extraction (ASE) in the Natural Products Industry.

The term natural products is a very broad classification and can be a part of Food and Beverage (Functional Foods) and Pharma/Life Science (Nutraceuticals) markets, and includes herbal health care and Traditional Chinese Medicine (TCM). Many scientists and organizations are working to understand what compounds are present in plant matrices that potentially cause the observed pharmacological effects. One of the challenges facing these scientists is the isolation of these compounds from the plant matrices. Often, traditional extraction methods produce extracts that are too complicated to analyze or characterize well. ASE® instruments increase the productivity of labs engaged in this field of research and produce extracts that are easier to analyze and characterize. Because water can be used as the extraction solvent, ASE instruments can also mimic one of the traditional modes of administration—seeping in hot water. Clearly, ASE technology is a good fit for any scientist or laboratory involved in isolating and studying natural products.

Features Unique to ASE	Benefits to Natural Product Extraction
True Walk-Away Automation	Overnight Extraction = Higher Productivity
Extract Same Sample with Different Solvents	Fractionation Extractions
Flow-Through Design	Removal of Interferences During Extraction
Acid/Base Extraction Capabilities	Hydrolyzation of Complex Matrices
Combined Static/Dynamic Extractions	Exhaustive Extraction = Increased Recoveries
Automatic Rinsing	System Rinsed with Any Solvent Automatically

## ASE Features and Benefits

### True Walk-Away Automation Details

This feature is very beneficial to labs extracting natural products because generally they have higher sample throughput, and preparation can be the bottleneck in the process. Using an ASE system, an entire rack of 24 can be set up at the beginning of the day and another rack of 24 samples can be set up to run overnight with extracts that are ready for analysis the first thing in the morning. With minimal user intervention, 48 samples, or more, per day can be extracted and ready for analysis. **No other sample preparation process available offers this type of automation!**

### Fractionation Extraction Details

Being able to extract the same sample into different collection vials allows the operator to use different solvents to extract different analytes from the same sample. For example, a plant sample can be extracted using weakly polar solvent first, removing the less polar compounds of interest, then the same sample can be extracted using a more strongly polar solvent to remove the more polar compounds of interest, with each extract collected in a separate vial for easy analysis. This makes the extracts easier to analyze because there are fewer compounds in each extract. **This feature is unique to ASE and cannot be duplicated by any other extraction method available!**

### Flow-Through Design Details

Because the ASE extraction cells are designed for the solvent to enter the top and exit the bottom of the cell with the analytes, the user can add various adsorbents to the bottom of the cell to remove unwanted co-extractables during the extraction process. This helps eliminate post extraction cleanup steps and accelerates the entire sample preparation process. This is very important to a laboratory performing extraction on natural products. Many times these samples are more difficult to extract and the extracts must be cleaned or filtered prior to analysis. With the ASE system, this labor-intensive step of the sample preparation process can be eliminated.

### Acid/Base Extraction Capabilities Details

Many times, natural product samples can be difficult to extract due to the complexity of the plant-based matrix. The acid- and base-resistant capabilities of the ASE system offer the customer some beneficial options. First, the samples can be hydrolyzed using a strong acid or base such as 8 M HCl or NaOH. This breaks down the plant matrix and allows the extraction solvent to reach the analytes of interest. The second option allows acidified or base-treated solvent to be pumped into the cell as the extraction is performed.

**Again, ASE is the only extraction method that offers this option!**

## Dynamic and Static Extraction Details

ASE is the only extraction method that provides both a dynamic and static extraction in the same run! Dynamic extraction is defined as the ability to introduce fresh solvent during the extraction process. This is important because it ensures that the extraction solvent will not become oversaturated with the analyte, decreasing its ability to remove more analyte. Static extraction is defined as holding the extraction solvent and sample for a set period of time to maximize the solubility of the analytes. Performing both dynamic and static extractions is what defines ASE as an exhaustive extraction technique and provides maximum analyte recoveries. This is important to laboratories extracting natural products to determine the quantity of active ingredients present in each sample.

## Automatic Rinsing Details

The automatic rinse function of the ASE system allows the user to set up different batches of samples for the same extraction run using different solvents for each batch. The system will automatically change solvents and rinse the entire system with the next solvent to be used, with no need for user intervention. This is ideal for laboratories extracting natural products, as many of the samples are different matrices and a run of 24 samples can contain any number of combinations of sample matrices and/or extraction methods.

ASE is the only extraction method to offer automatic rinsing!

# Application Notes

## AB 102: Determination of Aucubin, Geniposide, and Pinoselin Diglucoside in *Cortex Eucommiae* Using ASE and HPLC

**Overview:** This application brief (AB) describes an efficient method for the determination of aucubin, geniposide, and pinoselin diglucoside in *Cortex eucommiae* using Accelerated Solvent Extraction (ASE) and an UltiMate® 3000 HPLC system. The ASE method for the extraction of *C. eucommiae* replaced the soxhlet extraction method in the Chinese Pharmacopoeia Edition.

<http://www.dionex.com/en-us/webdocs/77703-AB102-ASE-LC-Cortex-18Aug09-LPN2192-01.pdf>

## AN 192: Rapid Analysis of Ginseng Using Accelerated Solvent Extraction and High Performance Liquid Chromatography

**Overview:** Here, the authors describe a method that combines ASE extraction with a 25 min HPLC separation for the analysis of 15 ginsenosides. This method is suitable for analyzing Asian ginseng, American ginseng, and notoginseng.

[http://www.dionex.com/en-us/webdocs/61830-AN192\\_ASE\\_HPLC\\_Ginseng\\_29Aug07\\_LPN1965.pdf](http://www.dionex.com/en-us/webdocs/61830-AN192_ASE_HPLC_Ginseng_29Aug07_LPN1965.pdf)

## AN 207: Chromatographic Fingerprinting of *Flos Chrysanthemi* Indici Using HPLC

**Overview:** In this application note, the authors establish an HPLC chromatographic fingerprint of *F. chrysanthemi* based on discriminating the characteristic peaks of chlorogenic acid and flavonoids (luteolin-7-o-glucoside, linarin, luteolin and apigenin; structures shown in Figure 1). An ultrasonic extraction technique is also compared to Accelerated Solvent Extraction for isolation of the target components from the samples.

<http://www.dionex.com/en-us/webdocs/71823-AN207-HPLC-Flos%20Chrysanthema-14April09-LPN2092.pdf>

## AN 232: Determination of Anthraquinones and Stilbenes in Giant Knotweed Rhizome by HPLC with UV Detection

**Overview:** An efficient HPLC method that determines the eight main active components of giant knotweed rhizome in a single injection (anthraglycoside A, anthraglycoside B, emodin, physcion, rhein, chrysophanol, resveratrol, and polydatin.) is described here. Samples were extracted using ultrasonic extraction and the ASE 200 Accelerated Solvent Extractor respectively, with satisfactory results obtained for both extraction methods.

<http://www.dionex.com/en-us/webdocs/77633-AN232-LC-Anthraquinones-Knotweed-21Aug2009-LPN2280-01.pdf>

## AN 335: Accelerated Solvent Extraction (ASE) of Active Ingredients from Natural Products

**Overview:** This application note describes the use of ASE for extraction of commercially available nutritional supplements; specifically hypericin from *Hypericum perforatum* (St. John's Wort) and berberine from *Hydrastis canadensis* (Goldenseal root). The current methods used for extraction of this product are Soxhlet and sonication, both of which are time- and solvent-intensive. Here, ASE extracts were generated in 14 minutes per sample using 15 mL of solvent and were immediately ready for processing and analysis.

<http://www.dionex.com/en-us/webdocs/4476-AN335.pdf>

## AN 346: Totally Automated Sample Preparation Using Accelerated Solvent Extraction (ASE) Coupled with the Gilson ASPEC®: The Determination of Dianthrone in St. John's Wort

**Overview:** Here, the authors report on the use of a coupled ASE-ASPEC system as an automated extraction and extract manipulation device for the determination of dianthrone in St. John's Wort (*Hypericum perforatum*). The results using the coupled system are compared to those obtained using a traditional sonication procedure followed by manual SPE cleanup.

[http://www.dionex.com/en-us/webdocs/4521-AN346\\_V13.pdf](http://www.dionex.com/en-us/webdocs/4521-AN346_V13.pdf)

## **AN 357: Extraction of Phenolic Acids from Plant Tissue Using Accelerated Solvent Extraction (ASE)**

**Overview:** The extraction of phenolic acids from two different plants (eggplants and black cohosh) using ASE is the focus of this method. ASE is able to extract phenolic acids from plant tissue more efficiently than traditional extraction techniques while saving time and solvent.

[http://www.dionex.com/en-us/webdocs/40398-AN357\\_V31\\_releasedJC120806.pdf](http://www.dionex.com/en-us/webdocs/40398-AN357_V31_releasedJC120806.pdf)

## **AN 362: Extraction of Herbal Marker Compounds Using Accelerated Solvent Extraction (ASE) Compared to Traditional Pharmacopoeia Protocols**

**Overview:** The use of the ASE system for solvent extraction of five marker compounds from herbal preparations followed by analysis by HPLC is reported here. The results are compared to traditional pharmacopoeia extraction methods.

<http://www.dionex.com/en-us/webdocs/69956-AN362-ASE-Marker-Compounds-Supplements-14Jan09-LPN2144.pdf>

## **Literature References**

4. F. Kawamura, Y. Kikuchi, T. Ohira, M. Yatagai.  
*ASE of paclitaxel and related compounds from the bark of taxus cuspidata*. Journal of Natural Products. 1999, 62, 244–247.
58. T-S. Peng, S. Cao, R. Davis, D. Johnston.  
*ASE of natural products*. 111th AOAC International Meeting. 1997:7–11 September, Paper 1102, .
59. B. Benthin, H. Danz, M. Hamburger.  
*Pressurized liquid extraction of medicinal plants*. Journal of Chromatography A. **1999**: 837, 211–219.
60. C.T. da Costa, S.A. Margolis, B.A. Benner, Jr., D. Horton.  
*Comparison of methods for extraction of flavanones and xanthenes from the root bark of the Osage orange tree using liquid chromatography*. Journal of Chromatography A. 1999: 831, 167–178.
139. D.F. Hein, H.E. Hummel.  
*Extraction of azadirachta indica ingredients: Comparison of soxhlet reflux extractor, tissue mixer, and accelerated solvent extraction*. Mededelingen-Faculteit Landbouwkundige en Toegepaste Biologische Wetenschappen (Universiteit Gent). 1999: 64,3a,185–195.
194. R.M. Alonso-Salces, E. Korta, A. Barranco, L.A. Berrueta , B. Gallo, F. Vincente.  
*Determination of polyphenolic profiles of basque cider apple varieties using ASE*. Journal of Agricultural Food Chemistry. **2001**: 49, 3761–3767.
249. E.S. Ong, S.M. Len.  
*Pressurized hot water extraction of berberrubine, baicalin and glycyrrhizin in medicinal plants*. Analytica Chimica Acta. **2003**: 482, 81–89.
289. A.L. Dawidowicz, D. Wianowska, J. Gawdzik, D.H. Smolarz.  
*Optimization of ASE conditions for the HPLC determination of rutin and isoquercitrin in sambucus nigra L*. Journal of Liquid Chromatography and Related Technologies. 2003:26,14, 2381–2397.
290. C.W. Huie.  
*A review of modern sample-preparation techniques for the extraction and analysis of medicinal plants*. Anal. Bioanal. Chem. **2002**, 23–30.
302. J.R. Denery, K. Dragull, C.S. Tang, Q.X. Li.  
*Pressurized fluid extraction of carotenoids from haematococcus pluvialis and dunaliella salina and kavalactones from piper methysticum*. Analytica Chimica Acta. **2003**, 175–181.
304. M. Palma, Z. Pineiro, C.G. Barroso.  
*Stability of phenolic compounds during extraction with superheated solvents*. Journal of Chromatography A. **2001**, 921, 169–174.
306. Z.Y. Ju, L.R. Howard.  
*Effects of solvent and temperature on pressurized liquid extraction of anthocyanins and total phenolics from dried red grape skin*. Journal of Agricultural and Food Chemistry. **2003**: 51, 5207–5213.
308. R. Wahlen, P. Norris, T. Catterick, M.D. Scrimshaw, J.N. Lester.  
*Comparison of microwave extraction and accelerated solvent extraction (ASE) for arsenic speciation in environmental samples and herbal medicine followed by HPLC/ICP-MS analysis*. LGC. **2002**.
314. Z. Pineiro, M. Palma, C.G. Barroso.  
*Determination of catechins by means of extraction with pressurized liquids*. Journal of Chromatography A. **2004**: 1026, 19–23.
411. F.Q. Yang, S.P. Li, Y. Chen, S.C. Lao, Y.T. Wang, T.T.X. Dong, K.W.K. Tsim.  
*Identification and quantitation of eleven sesquiterpenes in three species of curcuma rhizomes by pressurized liquid extraction and gas chromatography-mass spectrometry*. Journal of Pharmaceutical and Biomedical Analysis. **2009**: 39, 552–558.
416. M. Herrero, D. Arraez-Roman, A. Segura, E. Kenndler, B. Gius, M. Augusta Raggi, E. Ibanez, A. Cifuentes.  
*Pressurized liquid extraction-capillary electrophoresis-mass spectrometry for the analysis of polar antioxidants in rosemary extracts*. Journal of Chromatography A. **2005**: 1084, 54–62.
431. S. Mukhopadhyay, D.L. Luthria, R.J. Robbins.  
*Optimization of extraction process for phenolic acids from black cohosh (cimicifuga racemosa) by pressurized liquid extraction*. Journal of the Science of Food and Agriculture. **2006**: 86, 156–162.
432. D.L. Luthria, S. Mukhopadhyay.  
*Influence of sample preparation on assay of phenolic acids from eggplant*. Journal of Agricultural and Food Chemistry. **2006**: 54, 41–47.
435. A. Smelcerovic, V. Verma, M. Spiteller, S.M. Ahmad, S.C. Puri, G.N. Qazi.  
*Phytochemical analysis and genetic characterization of six hypericum species from Serbia*. Phytochemistry. **2006**: 67, 171–177.

436. V.F. Peres, J. Saffi, M.I.S. Melecchi, F.C. Abad, M.M. Martinez, E.C. Oliveira, R.A. Jacques, E.B. Caramao.  
**Optimization or pressurized liquid extraction of piper gaudi-chaudianum kunth leaves.** Journal of Chromatography A. **2006:** 1105, 148–153.
437. Z. Pineiro, M. Palma, C.G. Barroso.  
**Determination of trans-resveratrol in grapes by pressurized liquid extraction and fast high-performance liquid chromatography.** Journal of Chromatography A. **2006:** 1110, 61–65.
438. J. Zhao, S.P. Li, F.Q. Yang, P. Li, Y.T. Wang.  
**Simultaneous determination of saponins and fatty acids in ziziphus jujuba (suanzaoren) by high performance liquid chromatography-evaporative light scattering detection and pressurized liquid extraction.** Journal of Chromatography A. **2006:** 1108, 188–194.
439. Y. Jiang, S.P. Li, H.T. Chang, Y.T. Wang, P.F. Tu.  
**Pressurized liquid extraction followed by high-performance liquid chromatography for determination of seven active compounds in cortex dictamni.** Journal of Chromatography A. **2006:** 1108, 268–272.
452. H. Fan, S.P. Li, J.J. Xiang, C.M. Lai, F.Q. Yang, J.L. Gao, Y.T. Wang.  
**Qualitative and quantitative determination of nucleosides, bases and their analogues in natural and cultured Cordyceps by pressurized liquid extraction and high performance liquid chromatography-electrospray ionization tandem mass spectrometry (HPLC-ESI-M).** Analytica Chimica Acta. **2006:** 567, 218–228.
455. R. Anand, N. Verma, D.K. Gupta, S.C. Puri, V.K. Sharma, G.N. Qazi.  
**Comparison of extraction techniques for extraction of bioactive molecules from hypericum perforatum L. plant.** Journal of Chromatography A. **2005:** 43, 530–531.
456. S. Puri, G. Handa, V.K. Gupta, A.S. Shawl, O.P. Suri, G.N. Qazi.  
**Preparative high-performance liquid chromatographic separation of naphthodianthrone from St. John's wort.** Journal of Chromatographic Science. **2006:** 44, 177–180.
463. J.B. Wan, C.M. Lai, S.P. Li, M.Y. Lee, L.Y. Kong, Y.T. Wang.  
**Simultaneous determination of nine saponins from Panax notoginseng using HPLC and pressurized liquid extraction.** Journal of Pharmaceutical and Biomedical Analysis. **2006:** 41, 274–279.
470. A. Sae-Yun, C. Ovatlarnporn, A. Itharat, R. Wiwattanapatapee.  
**Extraction of rotenone from Derris elliptica and Derris malaccensis by pressurized liquid extraction compared with maceration.** Journal of Chromatography A. **2006:** 1125, 172–176.
485. S.C. Lao, S.P. Li, K.K.W. Kan, P. Li, J.B. Wan, Y.T. Wang, T.T.X. Dong, K.W.K. Tsim.  
**Identification and quantification of 13 components in Angelica sinensis (Danggui) by gas chromatography-mass spectrometry coupled with pressurized liquid extraction.** Analytica Chimica Acta. **2004:** 526, 131–137.
486. M. Palma, Z. Pineiro, C.G. Barroso.  
**In-line pressurized-fluid extraction-solid-phase extraction for determining phenolic compounds in grapes.** Journal of Chromatography A. **2002:** 968, 1–6.
488. P. Li, S.P. Li, S.C. Lao, C.M. Fu, K.K.W. Kan, Y.T. Wang.  
**Optimization of pressurized liquid extraction for Z-ligustilide, Z-butylidenephthalide and ferulic acid in Angelica sinensis.** Journal of Pharmaceutical and Biomedical Analysis. **2006:** 40, 1073–1079.
490. D. Luthria, S. Mukhopadhyay, A.L. Kwansa.  
**A systematic approach for extraction of phenolic compounds using parsley (Petroselinum crispum) flakes as a model substrate.** Journal of the Science of Food and Agriculture. **2006:** 86, 1350–1358.
491. L. Jaime, J.A. Mendiola, M. Herrero, C. Soler-Rivas, S. Santoyo, F.J. Senorans, A. Cifuentes, A. Ibanez.  
**Separation and characterization of antioxidants from Spirulina platensis microalga combining pressurized liquid extraction, TLC, and HPLC-DAD.** J. Sep. Sci. **2005:** 28, 2111–2119.
500. M. Herrero, E. Ibanez, J. Senorans, A. Cifuentes.  
**Pressurized liquid extracts from spirulina platensis microalga determination of their antioxidant activity and preliminary analysis by micellar electrokinetic chromatography.** Journal of Chromatography A. **2004:** 1047, 195–203.
503. M. Herrero, P.J. Martin-Alvarez, F. Javier Senorans, A. Cifuentes, E. Ibanez.  
**Optimization of accelerated solvent extraction of antioxidants from Spirulina platensis microalga.** Food Chemistry. **2005:** 93, 417–423.
507. E. Ibanez, M. Herrero, P.J. Martin-Alvarez, F.J. Senorans, G. Reglero, A. Cifuentes.  
**Accelerated solvent extraction: A new procedure to obtain functional ingredients from natural sources.** ACS Symposium Series 926. Modern Extraction Techniques Food and Agricultural Samples. **2006:** 65–78.
508. M. Waldeback, F.J. Senorans, A. Fridstrom, K.E. Markides.  
**Pressurized fluid extraction of squalene from olive biomass.** ACS Symposium Series 926. Modern Extraction Techniques Food and Agricultural Samples. **2006:** 96–106.
509. F.Q. Yang, Y.T. Wang, S.P. Li.  
**Simultaneous determination of 11 characteristic components in three species of Curcuma rhizomes using pressurized liquid extraction and high-performance liquid chromatography.** Journal of Chromatography A. **2006:** 1134, 226–231.
510. J.B. Wan, F.Q. Yang, S.P. Li, Y.T. Wang, X.M. Cui.  
**Chemical characteristics from different parts of Panax notoginseng using pressurized liquid extraction and HPLC-ELSD.** Journal of Pharmaceutical and Biomedical Analysis. **2006:** 41, 1596–1601.



525. D.L. Luthria.

**Application of green chemistry principles for extraction of phytolipids and phenolic compounds.** Indian Journal of Chemistry. **2006:** 45B, 2291–2296.

527. M. Herrero, E. Ibanez, A. Cifuentes, g. Reglero, S. Santoyo.

**Dunaliella salina microalga pressurized liquid extracts as potential antimicrobials.** Journal of Food Protection. **2006:** 69,10, 2471–2477.

532. N.Y. Qin, F.Q. Yang, Y.T. Wang, S.P. Li.

**Quantitative determination of eight components in rhizome (Jianghuang) and tuberous root (Yujin) of Curcuma longa using pressurized liquid extraction and gas chromatography-mass spectrometry.** Journal of Pharmaceutical and Biomedical Analysis. **2007:** 43, 486–492.

543. J. Zhao, X-Q. Zhang, S-P Li, F-Q Yang, Y-T Wang, W-C Ye.

**Quality evaluation of Ganoderma through simultaneous determination of nine triterpenes and sterols using pressurized liquid extraction and high performance liquid chromatography.** J. Sep. Sci. **2006:** 29, 2609–2615.

546. Xu Xiao-Qin, Li Qing-Ling, Yuan Ji-Duan, Wang Shu-Gui, Wang Wen-Shen.

**Determination of three kinds of chloroacetanilide herbicides in Radix Pseudostellariae by accelerated solvent extraction and gas chromatography-Mass Spectrometry.** Chinese Journal of Analytical Chemistry. **2007:** 35,2, 206–210.

550. Chunhui Deng, Ning Liu, Mingxia Gao, Xiangmin Zhang.

**Recent developments in sample preparation techniques for chromatography analysis of traditional Chinese medicines.** Journal of Chromatography A. **2007:** 1153, 90–96.

563. Emma Warburton, Paul L Norris, Heidi Goenaga-Infante.

**Comparison of the capabilities of accelerated techniques for the quantification of kavalactones in piper methysticum (kava) roots by high performance liquid chromatography with ultra violet detection.** Phytochemical Analysis. **2007:** 18, 98–102.

564. Tobias Mohn, Olivier Potterat, Matthias Hamburger.

**Quantification of Active principles and pigments in leaf extracts of isatis tinctoria by HPLC/UV/MS.** Institute of Pharmaceutical Biology. **2007:**73, 151–156.

## Customer References

Names of scientists using ASE for natural products can be obtained by contacting the Salt Lake Technical Center directly or via email at [asesupport@dionex.com](mailto:asesupport@dionex.com).

ASE and UltiMate are registered trademarks of Dionex Corporation. ASPEC is a trademark of Gilson Incorporated. All other trademarks and registered trademarks are the property of Dionex Corporation.

Passion. Power. Productivity.



#### Dionex Corporation

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

#### North America

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

#### South America

Brazil (55) 11 3731 5140

#### Europe

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

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



LPN 2573 PDF 07/10  
©2010 Dionex Corporation