

Instruction of the Clean-up Process using *B-TeZ IAC Folic Acid 3ml*



Fast and Accurate Content Determination of Folic Acid in Vitamin Tablets, Liquid Vitamin Preparations, Cell Culture Extracts etc. by Combination of Immunoaffinity Chromatography and HPLC-UV

Principle:

Many methods of Folic Acid determination based on HPLC-UV (High Performance Liquid Chromatography-Ultraviolet) or HPLC-FLD (High Performance Liquid Chromatography with Postcolumn Fluorescence Derivatization) detection show low selectivity if problematic matrices are applied.

This method of determination of Folic Acid combines the high selectivity of immunoaffinity columns (compared to other separation materials (e.g. SAX)) with its potential to concentrate elute through use of purification by a HPLC column.

Sample Preparation:

Folic Acid samples are to be extracted and analyzed with the method of E.S. Osseyi et al. [E.S. Osseyi, R.L. Wehling, J.A. Albrecht *J. Chromatogr. A* 1998; 826:235-240], e.g. vitamin tablets, liquid vitamin preparations, cell culture extracts. Example: 2g vitamin containing tablets are dissolved in 50ml PBS. The resulting extract may be filtered through a 0.45µm membrane filter.

Enrichment Step IAC:

10ml extract (containing the quantity of folic acid from a 0.4g sample if above-mentioned sample preparation is followed) is diluted with a total volume of 30ml PBS and then applied in a reservoir on top of the *B-TeZ IAC Folic Acid 3ml* column. The optimal flow rate through the gel is between 1 to 3 ml/min.

Wash:

After whole sample has passed through the gel the latter is washed with 5ml of PBS. Remaining liquids in the gel are removed by applying either pressure from top of the column or pressure from the bottom.

Elution:

The sample reservoir on top of the *B-TeZ IAC Folic Acid 3ml* column is removed, and an appropriate vial is placed below the affinity column. The bounded folic acid is eluted by using a total volume of 3ml of methanol-phosphoric acid solution (99.8/0.2 v/v). The elution process

is performed in two steps. First, an amount of 1ml methanol is applied. Once this amount has passed through the column, there should be a waiting time of 30 seconds. After that, the second portion of 2ml of methanol is eluted through the column. The flow rate should lie below 3ml/min. The remaining methanolic solutions should be eluted by application of slight under- or overpressure. All methanolic fractions are unified to give the column elute.

The column elute may be injected into the HPLC directly or, if concentrations are very low, concentrated by evaporation at 50°C for 1h (e.g. using VLM evaporator), re-dissolved in HPLC solvent and finally injected into the system. For the latter case, please see the sample calculation in which the sample concentrate is re-dissolved in 1.0ml HPLC solvent.

Analytical Method:

- Machine: Shimadzu
- Column: Trentec Reprisil-Pur RP C18 120 ODS3 5µm 125x3,0mm with guard column
- Mobile Phase B: 0.03M potassium phosphate, pH 2.2-methanol (80/20 v/v)
- Gradient: 0.01min B 100%; 25min B 100% (isocratic)
- Flow Rate: 0.5ml/min
- Time of Analysis: 25min
- Injector Volume: 100µl
- Detection: λ ABS [nm]: 280nm.

Characteristics:

The linear range is 20ng to 500ng Folic Acid, per injection ($R^2=0.9946$). The limit of detection is 3 ng of folic acid per injection (3 times of signal/noise ratio). If the given dilution steps are followed, the folic acid contents of 0.5 to 12.5µg/g lie within the linear working range of the method. The lower limit of detection is 75ng/g of folic acid in the sample. The lower limit of detection may be reduced by enhancement of folic acid in the sample.

Recovery rates are >85% when folic acid in buffer mixtures is analysed in the range of 0.1 to 5µg per IAC.

Instruction of the Clean-up Process using *B-TeZ IAC Folic Acid 3ml*



Example Sample Calculation:

$\frac{2\text{g Sample}}{50\text{ml Extraction Solvent}} \times \frac{10\text{ml Extract}}{1\text{ml}} \times 0.1\text{ml injector volume} = 0.04\text{g Sample Equivalents}$

$\frac{\# \mu\text{g injected Folic Acid}}{\text{Sample Equivalents [g]}} = \mu\text{g/g Folic Acid}$

Buffer and Chemicals:

Acidic Elution Solvent:

methanol-phosphoric acid (99.8/0.2 v/v)

Dissolve 200µl phosphoric acid (95%) in 100ml methanol

HPLC-Solvent

0.03M potassium phosphate, pH 2.2-methanol (80/20 v/v)

Dissolve 4.1g KH₂PO₄ in 800ml deionized water. Adjust to pH 2.2 with phosphoric acid (95%). Add 200ml methanol. Degass with helium.

Chemicals:

- methanol, HPLC grade
- deionized water
- dipotassium hydrogenphosphate, >98%
- potassium dihydrogenphosphate, >98%
- sodium chloride
- phosphoric acid, 95%

Consumables:

- *B-TeZ IAC Folic Acid 3ml* column [BTFS319005]

Standard:

- Folic Acid (98%) [Sigma F-7876]

Evaporation:

- nitrogen gas 5.0 [Air Liquide M55763810] (to evaporate IAC-eluate)

Apparatus:

- HPLC; Shimadzu; pump: LC-6A (2 pieces); autosampler: SIL 6B; fluorescence detector: RF-10AXL; data handling: CLASS LC10

- Vakuum SPE Manifold (BAKER spe-24G Column Processor – process up to 24 samples) [J.T. Baker 7208]

- Evaporator (with tripod) [VLM EVA EC1-S]

Instruction of the Clean-up Process using **B-TeZ IAC Folic Acid 3ml**



Manufactured by:

BioTez Berlin-Buch GmbH
Robert-Rössle-Str. 10, 13125 Berlin Germany
Phone +49 30 9489-2130 • Fax +49 30 9494-509
e-mail info@biotez.de • www.biotez.de

distributed in the US/Canada by:



EAGLE BIOSCIENCES, INC.
20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 • FAX: 617-419-1110
WWW.EAGLEBIO.COM • info@eaglebio.com

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.