

# Instruction of the Clean-up Process using *B-TeZ IAC Biotin 3ml*



## Fast and Accurate Content Determination of Biotin in Vitamin Tablets, Liquid Vitamin Preparations, Cell Culture Extracts, etc. by Combination of Affinity Chromatography, HPLC-FLD and Post Column Derivatisation with Fluorescein-Streptavidin Conjugate as Fluorophor

### Principle:

Many methods of biotin determination based on HPLC-UV (High Performance Liquid Chromatography Ultraviolet) detection may show either low sensitivity or low selectivity. This depends on the dilution factor of the matrices and if problematic matrices are applied.

This method of content determination of Biotin combines the high selectivity of affinity columns with its potential to concentrate elute and use high sensitivity biotin detection by post column labelling with fluorescein-streptavidin conjugate.

### Sample Preparation:

Biotin samples are to be extracted and analyzed with the method of Bachas et al. [N.G. Hentz, L.G. Bachas Methods Enzymol. 1997; 279:275-86], e.g. vitamin tablets, liquid vitamin preparations, cell culture extracts. Example: 25g vitamin containing tablets are dissolved in 100ml PBS. The resulting extract may be filtered through a 0.45µm membrane filter.

### Enrichment Step IAC:

4ml extract (containing the quantity of Biotin from a 1g sample of above-mentioned sample preparation is followed) **is diluted with a total volume of 20ml PBS** and then applied in a reservoir on top of the BioTeZ-Immunoaffinity Column. **The optimal flow rate through the gel is between 1 to 3ml/min.**

### Wash:

After the 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 BioTeZ-Immunoaffinity Column is removed, and an appropriate vial is placed below the affinity column. The bounded biotin is **eluted by using a total volume of 3ml of HPLC grade methanol**.

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 are 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 0.4ml HPLC solvent.

### Analytical Method:

- Machine: Shimadzu
- Column: Trentec Reprosil-Pur RP C18 120 ODS3 5µm; 125x3,0mm with guard column
- Mobile Phase A: methanol /water (85:15 v/v) (use only for cleaning purposes at the beginning and at the end of analytical series)
- Mobile Phase B: 0.1M potassium phosphate, pH 7.0-methanol (85/15 v/v)
- Gradient: 0.01min B 100%; 12min B 100%
- Flow Rate: 0.4ml/min
- Time of Analysis: 12min
- Injector Volume: 100µl
- Detection:  $\lambda_{EX}$  [nm]: 495nm;  $\lambda_{EM}$  [nm]: 518nm.
- Post column derivatization: A solution of Fluorescein (FITC)-Streptavidin-Conjugate (e.g. BioTeZ-FITC-Streptavidin) of protein concentration of 2µg/ml in 0.1M potassium phosphate, pH=8.2; flow rate: 0.2ml/min.

# Instruction of the Clean-up Process using *B-TeZ IAC Biotin 3ml*



## Characteristics:

The HPLC measuring range is linear of 5ng to 40ng Biotin per injection ( $R^2=0.994$ ). The limit of detection is 2ng of biotin per injection (three times of signal/noise ratio). If the given dilution steps are obeyed, the Biotin contents of 20 to 160 ng/g lie within the linear working range of the method. If the contents of used samples are higher than the cited range, extracts or the IAC column elutes should be diluted in a suitable manner. The lower limit of quantization is 10ng/g of Biotin in the sample.

**Recovery rates** are >85% when Biotin contents in buffer mixtures are analysed in the range of 0.1 to 5µg per IAC.

## Example Sample Calculation:

$$\frac{25\text{g Sample}}{100\text{ml Extraction Solvent}} \times \frac{4\text{ml Extract}}{0.4\text{ml}} \times \frac{0.1\text{ml}}{\text{injector volume}} = 0.25\text{g Sample Equivalents}$$

$$\frac{\text{\# } \mu\text{g injected Biotin}}{\text{Sample Equivalents [g]}} = \mu\text{g/g Biotin in e.g. multivitamin tablets}$$

## Buffer and Chemicals:

### **Phosphate Buffered Saline pH 7.4 (= PBS)**

1.24g  $\text{KH}_2\text{PO}_4$   
7.27g  $\text{K}_2\text{HPO}_4$   
8.76g NaCl

Dissolve in 1L deionized water. If necessary adjust pH to 7.4

### **HPLC-Solvent**

0.1M potassium phosphate, pH 7.0-methanol (85/15 v/v)

Dissolve 13.6g  $\text{KH}_2\text{PO}_4$  in 850ml deionized water. Adjust to pH 7.0 with 1M NaOH. Add 150ml methanol. Degas with helium.

0.1M potassium phosphate, pH 8.2

Take 200ml 0.1M potassium phosphate, pH 7.0. Adjust to pH 8.2 with 1M NaOH. Degas with helium.

methanol /water (85:15 v/v)  
(HPLC Column Cleaning)

Mix 85ml methanol and 15ml deionized water. Degas with helium.

### **Chemicals:**

- acetonitrile, HPLC grade
- methanol, HPLC grade
- deionised water
- dipotassium hydrogenphosphate, >98%
- potassium dihydrogenphosphate, >98%
- sodium chloride

### **Consumables:**

- B-TeZ Biotin Affinity Column

### **Standard:**

- D(+)-Biotin, 99% [SigmaAldrich B4501]

# Instruction of the Clean-up Process using **B-TeZ IAC Biotin 3ml**



## Flourescence Label for Biotin:

- BioTeZ Streptavidin-FITC-Conjugate  
(Order Code: StrAv-FITC)

## Evaporation:

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

## Apparatus:

- HPLC; Shimadzu; pump: LC-6A (2 pieces); auto sampler: SIL 6B; fluorescence detector: RF-10AXL; data handling: CLASS LC10
- Vacuum SPE Manifold (BAKER spe-24G Column Processor – process up to 24 samples [J.T. Baker 7208])
- Evaporator (with tripod) [VLM EVA EC1-S]



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](mailto:info@biotez.de) • [www.biotez.de](http://www.biotez.de)

distributed in the US/Canada by:



**EAGLE**  
BIOSCIENCES

### EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063  
Phone: 617-419-2019 • FAX: 617-419-1110  
[WWW.EAGLEBIO.COM](http://WWW.EAGLEBIO.COM) • [info@eaglebio.com](mailto: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](mailto:info@eaglebio.com) or at 866-411-8023.*