

Instruction of the Clean-up Process Using *B-TeZ IAC Aflatoxin M₁ 3ml*



Clean-up of Milk of Milk Products containing Aflatoxin M₁ (AFM₁) Via Immunoaffinity Chromatography and Subsequent HPLC Analysis with Postcolumn Derivatization with Perbromide (for joint determination of Aflatoxin B₁)

Principle:

Immunoaffinity chromatography stands for an efficient enrichment process of aflatoxins with concomitant effective separation from matrix components. The eluate is best suited for determination by HPLC. By varying washing process, that means exchange PBS buffer with 0.2M ammonium acetate solution, GC can also be applied.

Sample Preparation of Milk:

According to the method of Shundo et al.¹ the milk to be tested is first centrifuged by 1540g for 15 minutes. The cream is then separated using e.g. a separation funnel. If separation is incomplete, the milk sample should be filtered to avoid that remaining fatty particles block the column in next step below.

Enrichment Step IAC:

50ml skimmed milk are diluted with 5ml 10X PBS (= 10fold concentrated 50mM PBS buffer) and then applied in a reservoir on top of the *B-TeZ IAC Aflatoxin M₁ 3ml* column. Rate of flow through the affinity gel is 1 to 3 ml/min.

Caution! Be aware that no big air bubbles are neither in the gel nor between gel and luer lock outlet of column which prevent a permanent flow or necessary exchange of matter.

Depending on application and on expected contents, larger or smaller extract aliquots can be applied. In such cases the sample calculation (see below) must be adapted.

If samples are to be prepared simultaneously, manifold of J.T. Baker for 12 samples has proven of value.

Wash:

After whole sample has passed through the gel the latter is washed with 5ml of 50mM PBS. Remaining liquids in the gel are removed by applying either pressure from top of the column or underpressure from bottom.

Elution:

Sample reservoir on top of the *B-TeZ IAC Aflatoxin M₁ 3ml* column is removed and a appropriate vial is placed below the affinity column. The bound toxins are eluted by using a total of 1.5ml of methanol as elution solvent. The elution process is performed in two steps to ensure complete release of analytes. First, a volume of 0.5ml elution solvent is applied. After that volume has passed through column half a minute is waited before the second portion of 1ml of elutions solvent is eluted through the column. Remaining solvent solutions should be eluted by application of slight under- or overpressure. All methanolic fractions are unified to give the column eluate.

The column elute may be injected into the HPLC directly or in case concentrations are low it may be concentrated by evaporation (e.g. using VLM evaporator at 50°C under a permanent stream of nitrogen).

Caution: If you follow concentration by evaporation, add 100µl of acetic acid/water (50/50 v/v) as keeper. To avoid loss of analytes, stopp concentration at a small volume of residue, e.g. 50 to 100µl.

Instruction of the Clean-up Process Using **B-TeZ IAC Aflatoxin M₁ 3ml**



The residue then is redissolved in HPLC solvent, e.g. in 400µl in example calculation see below, and a aliquot is finally injected into the system.

IAC Column Characteristics:

A) Working Range and Recovery Rates of B-TeZ IAC Aflatoxin M₁ 3ml Column:

Working Range of Column:	0.04 – 50ng AFM ₁ per IAC
Zero Contamination of Column:	<0.04ng (LOD of HPLC-FLD method)
Guaranteed Recovery Rates for AFM ₁ within the Working Range ^(*) :	>85%

^(*) Recovery rates are confined to solvent content of diluted extract below 20% methanol or 10% acetonitrile.

B) Cross Reactivities^() of B-TeZ IAC Aflatoxin M₁ 3ml Column:**

Aflatoxin M ₁ (AFM ₁):	100%
Aflatoxin B ₁ (AFB ₁):	98%
Aflatoxin B ₂ (AFB ₂):	99%
Aflatoxin G ₁ (AFG ₁):	50%
Aflatoxin G ₂ (AFG ₂):	98%

^(**) Ratio of IAC recovery rates if a quantity of 5ng Aflatoxin total per column is analyzed

C) Capacity^(*) of B-TeZ IAC Aflatoxin M₁ 3ml Column:**

Maximum Column Capacity:	0.9µg AFM ₁
--------------------------	------------------------

^(***) An excess of AFM₁, e.g. 3µg, in a small volume of 2ml PBS is incubated with the IAC for 5 minutes; then the IAC is washed with 2ml PBS and the nonbonded fraction is analyzed. The difference of added analyt and nonbonded analyt equals maximum column capacity.

Analytical Method:

HPLC: Shimadzu; Column: Trentec Reprosil-Pur RP C18 120 ODS3 5µm; 125x3,0mm with guard column; Mobile Phase A: methanol / deionized water (85/15 v/v); Mobile Phase B: methanol / acetonitrile / deionized water (18/18/64 v/v/v); Gradient: 0,01 min B 100 %; 16 min B 100 %; 17 min B 0 %; 19 min B 0 %; 20 min B 100 %; Flow Rate: 0.5ml/min; Time of Analysis: 30min; Injector Volume: 100µl

Post Column Derivatization (only advisable in case of concomitant aflatoxin B₁ analysis): 32 ppm pyridinium hydrobromide perbromide in dioxan/ deionized water (0.1/99.9, v/v); Flow Rate: 0.2ml/min; reaction capillary put between end of column and detector by T-device: PEEK-capillary, 1/16'' x 0.25 mm ID; length: 40 cm

Fluorescence-Detection: λ_{EX} [nm]: 362nm; λ_{EM} [nm]: 440nm.

Temperature: Machine and eluents are at room temperature. Eluents are degassed with helium gas.

Example Sample Calculation of AFM₁ content:

(Calculation of AFB₁ is analogous)

A) Calculation of Sample Gramm Equivalents per HPLC injection:

$$\frac{50\text{g Sample}}{55\text{ml Extraction Mixture}} \times \frac{55\text{ml Extract}}{0.4\text{ml}} \times \frac{0.1\text{ml}}{\text{injector volume}} = 12.5\text{g Sample Equivalents}$$

Instruction of the Clean-up Process Using *B-TeZ IAC Aflatoxin M₁ 3ml*



B) Calculation of Aflatoxin M₁ contamination of examined commodity in ng/g:

$\frac{\text{\# ng injected AFM}_1}{\text{Sample Equivalents [g]}} = \text{ng/g AFM}_1 \text{ in commodity}$
--

Buffer, Chemicals, Apparatus and Literature:

10X PBS buffer:

12.4g KH₂PO₄
72.7g K₂HPO₄
87.6g NaCl

Dissolve in 1L deionised water. If necessary
adjust pH to 7.2 (± 0.2) with 1N NaOH or 1N HCl

Phosphate Buffered Saline pH 7.4 (= 50mM PBS):

Dilute 100ml 10X PBS with deionized water to a
final volumen of 1L. Control pH to 7.4 (± 0.2)

Chemicals:

- acetonitrile, HPLC grade
- methanol, HPLC grade
- acetic acid, 100% ultrapure

Consumerables:

- B-TeZ IAC Aflatoxin M₁ 3ml*
[BTAFM311005]

Elution Solvent:

Methanol

- deionized water

- dipotassium hydrogenphosphate, >98%
- potassium dihydrogenphosphate, >98%
- sodium chloride

Mycotoxin Standard:

- Aflatoxin M1 [e.g. Biopure]

Evaporation:

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

Keeper:

acetic acid/water (50/50 v/v):
Mix 10ml acetic acid and 10ml deionized water

Postcolumn Derivatization:

- pyridine hydrobromide perbromide, >95%.
- dioxane, 99.5%

Reagent:

32 ppm pyridine hydrobromide perbromide in
dioxan / deionized water (0.1/99.9, v/v):

Partially dissolve 32mg pyridine hydrobromide
perbromide in 1ml dioxane and pipet the oily
suspended liquid into 1L degassed deionized
water. Mix thoroughly.

Instruction of the Clean-up Process Using ***B-TeZ IAC Aflatoxin M₁ 3ml***



Apparatus:

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

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

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

BioTeZ Berlin-Buch GmbH

www.biotetz.de

¹ "AFLATOXIN M1 IN MILK BY IMMUNOAFFINITY COLUMN CLEANUP WITH TLC/HPLC DETERMINATION" Luzia Shundo; Myrna Sabino, *Brazilian Journal of Microbiology* **2006**, 37,164-167



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

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.