

Instruction of the Clean-up Process Using **B-TeZ IAC T2/HT2 3ml**



Clean-up of Commodity Extracts of Food and Feed Samples containing T2/HT2 via Immunoaffinity Chromatography and Subsequent Determination by HPLC with Precolumn Derivatization using 1-Anthroylnitrile and 4-Dimethylaminopyridin

Principle:

This instruction of T2/HT2 determination in food and feed focuses on the enrichment step of extract using immunoaffinity column (IAC) and quantification with HPLC.

Accepted laboratory extraction methods could be maintained. Full performance of the IAC column is given if pronounced criteria regarding organic solvent tolerance, elution process of analyte and working range of column is followed.

Many pretreatment methods for T2/HT2 in food and feed show low sensitivity because of interfering substances if problematic matrices are applied.

This method of content determination of T2/HT2 combines the high selectivity of an immunoaffinity column (IAC) with its potential to concentrate elute and additional step of purification of labelled T2/HT2 by HPLC column.

As said before, this instruction focuses on the handling with the IAC column. For the commodity extraction step a literature method is given. Please see below. The given apparatus (e.g. HPLC system) might serve as example among other possibilities. For your convenience, an example HPLC method for the analysis of T2-toxin is given below.

Sample Preparation (Literature method given):

Samples which content of T2/HT2 are to be analyzed, e.g. cereals, are extracted by the method of Visconti et al.¹ using methanol-water (90/10 v/v) as extraction solvent. E.g. to 50g of sample are added a volume of 100ml of the extraction solvent and processed as cited.

Enrichment Step IAC:

2ml extract (1g sample equivalent if above mentioned example extraction is followed) is diluted with 18ml 10mM PBS, pH=7.2 resulting a total volume of 20ml and then applied in a reservoir on top of the **B-TeZ IAC T2/HT2 3ml** column.

If the example is followed the resulting organic solvent concentration of the diluted extract is 9% of methanol which is tolerated by the column. The flow rate is adjusted to lie between 1 to 3ml/min.

According to application and contents expected the applied extract volumes could vary. In case of very low contents even extract volumes of 20ml may be diluted with buffer and applied without significant loss of analyt as long as resulting pH is fairly neutral and alcohol or acetonitrile content lies under 15%.

If latter is not the case the extract must further be diluted with PBS until forementioned maximum allowed organic solvent content during enrichment step using the **B-TeZ IAC T2/HT2 3ml** column is reached.

Thus, to guarantee maximum efficiency in terms of maximum recovery rates using **B-TeZ IAC T2/HT2 3ml** column, please ensure that

1. organic solvent content in diluted extract which is applied on top of the column is not higher than 15%,
2. T2/HT2 load per column does not exceed a total of 1µg (see working range below).

Instruction of the Clean-up Process Using **B-TeZ IAC T2/HT2 3ml**



Wash:

After whole sample has passed through the gel the latter is washed with 5ml of 10mM 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 using **B-TeZ IAC T2/HT2 3ml** column is removed and an 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 methanol fractions are unified to give the column eluate.

The column eluate is concentrated to complete dryness (e.g. using VLM evaporater at 50°C under a permanent stream of nitrogen). The residue is treated with a mixture of 1-Anthroylnitrile and 4-Dimethylaminopyridin as reported¹ to give fluorescent labelled derivatives of T2 and HT2 which are then injected to the HPLC apparatus.

Column Characteristics:

A) Working Range and Recovery Rates of B-TeZ IAC T2/HT2 3ml Column:

Working Range of Column:	upto 1µg T2/HT2 per IAC
Zero Contamination of Column:	<1ng T2/HT2 (LOD of HPLC method)
Guaranteed Recovery Rates within the Working Range ^(*) :	
T2-Toxin:	>85%
HT2:	>85%

^(*) Recovery rates are confined to solvent content of diluted extract below 15% methanol or acetonitrile (see details under Enrichment Step).

B) Cross Reactivities^() of B-TeZ IAC T2/HT2 3ml Column:**

T2-Toxin:	85%
HT2:	100%
DON:	<1%

^(**) Recovery rates if a total of 1.5µg T2/HT2/DON (1:1:1 molar ratio) is analysed per column.

C) Capacity^(*) of B-TeZ IAC T2/HT2 3ml Column:**

Maximum Column Capacity:	3µg T2 Toxin
--------------------------	--------------

^(***) An excess of T2 Toxin, e.g. 5µg T2, 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 analyte and nonbonded analyte equals maximum column capacity.

Analytical Method:

HPLC: Shimadzu; **Column:** Trentec Select RP C8 120 ODS3 5µm; 125x3,0mm with guard column; **Mobile Phase A:** acetonitrile / water (90/30, v/v); **Mobile Phase B:** acetonitrile / water (60/40, v/v); **Gradient:** 0.01min B 70%; 8min B 70%; 8.1min B 0%; 20min B 0%; 20.1min B 70%; **Flow Rate:** 0.5ml/min; **Time of Analysis:** 30min; **Injector Volume:** 100µl; **Detection:** λ_{EX} [nm]: 370nm; λ_{EM} [nm]: 470nm.

Instruction of the Clean-up Process Using *B-TeZ IAC T2/HT2 3ml*



Overall Method Characteristics (IAC and HPLC):

Measuring range is linear of 5ng to 100ng T2 pro injection ($R^2=0.999$). Limit of detection is 1ng of T2 pro injection (3times of signal/noise ratio). HT2 characteristics are quite similar. If given dilution steps are obeyed T2/HT2 commodity contents totalling **0.05 to 1µg/g** lie within the linear working range of the method (included working range of IAC column). If contents of used samples are higher than cited upper range, smaller portion of extract should be applied on the IAC column and analysed to ensure maximum recovery rates.

Example Sample Calculation of T2 content:

(Calculation of HT2 content is analogous)

50g Sample -----	x	2ml Extract -----	x	0.1ml injector volume	= 0.1g Sample Equivalents
100ml Extraction Solvent		1ml			

# µg injected T2 -----	=	µg/g T2 in e.g. cereal
Sample Equivalents [g]		

Buffer, Chemicals, Apparatus and Literature:

Phosphate Buffered Saline pH 7.4 (= 10mM PBS)

0.25g KH_2PO_4
1.45g K_2HPO_4
8.76g NaCl

Dissolve in 1L deionized water. If necessary
adjust pH to 7.4 (± 0.3) with 1N NaOH or 1N HCl

Chemicals:

- acetonitrile, HPLC grade
- methanol, HPLC grade
- deionized water
- dipotassium hydrogenphosphate, >98%
- potassium dihydrogenphosphate, >98%
- sodium chloride
- nitrogen gas 5.0 [Air Liquide M55763810] (to evaporate IAC-eluate)

Consumables:

- ***B-TeZ IAC T2/HT2 3ml*** [Prod. No. BTTE316005]

Elution Solvent:

methanol

OH Derivatization:

- 1-Anthroyl Nitrile [WAKO 017-12101]

Dilution Solvent:

acetonitrile-water (40/60 v/v):

Instruction of the Clean-up Process Using **B-TeZ IAC T2/HT2 3ml**



- 4-Dimethylamino-pyridin [SIGMA-ALDRICH 29224-10G] Mix 40ml acetonitrile and 60ml water.
- Toluene [ROTH 7341.1]

Apparatus:

HPLC; Shimadzu; pump: LC-6A (2 pieces); autosampler: 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.biotez.de

¹ "Analysis of T-2 and HT-2 toxins in cereal grains by immunoaffinity clean-up and liquid chromatography with fluorescence detection", Angelo Visconti, Veronica Maria Teresa Lattanzio, Michelangelo Pascale, Miriam Haidukowski, *Journal of Chromatography A*, 1075 (2005) 151–158

Biodiagnostika
Oligonucleotide
Markierte Biomoleküle



BIOTEZ

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



EAGLE
BIOSCIENCES

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.