

Total Aflatoxin ELISA Kit ab285282

1 图像

概述

产品名称	Total Aflatoxin ELISA试剂盒
检测方法	Colorimetric
检测类型	Competitive
灵敏度	0.02 ng/ml
实验步骤	Multiple steps standard assay
产品概述	<p>This Total Aflatoxin ELISA Kit (ab285282, E4747) is based on the Competitive ELISA method. It can detect Aflatoxin (AF) in samples, such as grain, peanut, formula feed, etc.</p> <p>The microplate provided in this kit has been pre-coated with coupled antigen. During the reaction, AF in the samples or standard competes with coupled antigen on the solid phase supporter for sites of anti-AF antibody. Then Horseradish Peroxidase (HRP) conjugate is added to each micro plate well, and TMB substrate is added for color development.</p> <p>There is a negative correlation between the OD value of samples and the concentration of AF. The concentration of AF in the samples can be calculated by comparing the OD of the samples to the standard curve.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called E4747 Total Aflatoxin ELISA Kit. E4747-100 is the same size as the 96 test size of ab285282.</p>
平台	Pre-coated microplate (12 x 8 well strips)

性能

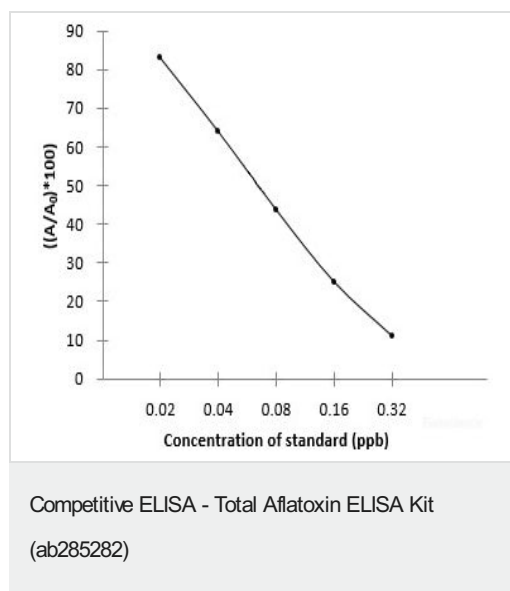
存放说明Store at +4°C. Please refer to protocols.

组件	96 tests
20X Wash buffer	1 x 40ml
Antibody Working Solution	1 x 5.5ml
ELISA Microplate	1 unit

组件	96 tests
HRP Conjugate	1 x 5.5ml
Plate Sealer	3 units
Standards (S1 - S6)	6 x 1ml
Stop Solution	1 x 6ml
Substrate Reagent A	1 x 6ml
Substrate Reagent B	1 x 6ml

相关性 Aflatoxins are naturally occurring mycotoxins that are produced by many species of *Aspergillus*.

图片



Example of standard curve using ab285282. This standard curve is for demonstrative purposes only. A standard curve must be run with each assay.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors