

Aflatoxin B1 (AFB1) ELISA Kit ab302933

概述

|      |                                                                                                                                                                                                                                   |   |      |              |
|------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|---|------|--------------|
| 产品名称 | Aflatoxin B1 (AFB1) ELISA试剂盒                                                                                                                                                                                                      |   |      |              |
| 检测方法 | Colorimetric                                                                                                                                                                                                                      |   |      |              |
| 精确度  | 批次内                                                                                                                                                                                                                               |   |      |              |
|      | 样品                                                                                                                                                                                                                                | n | Mean | SD           |
|      | Overall                                                                                                                                                                                                                           |   |      | CV%<br>< 10% |
|      | 批次间                                                                                                                                                                                                                               |   |      |              |
|      | 样品                                                                                                                                                                                                                                | n | Mean | SD           |
|      | Overall                                                                                                                                                                                                                           |   |      | CV%<br>< 10% |
| 样品类型 | Food samples                                                                                                                                                                                                                      |   |      |              |
| 检测类型 | Competitive                                                                                                                                                                                                                       |   |      |              |
| 灵敏度  | = 0.15 ng/ml                                                                                                                                                                                                                      |   |      |              |
| 范围   | 0.15 ng/ml - 4.05 ng/ml                                                                                                                                                                                                           |   |      |              |
| 检测时间 | 0h 25m                                                                                                                                                                                                                            |   |      |              |
| 实验步骤 | Multiple steps standard assay                                                                                                                                                                                                     |   |      |              |
| 产品概述 | Aflatoxin B1 (AFB1) ELISA Kit (AB302933), previously known as Aflatoxin B1 (AFB1) ELISA Kit (Catalog # K4208), is a competitive ELISA kit for the quantitative measurement of Aflatoxin B1 in feedstuff, feed, grain, edible oil. |   |      |              |
|      | Detection limit:                                                                                                                                                                                                                  |   |      |              |
|      | Edible oil 1 µg/kg                                                                                                                                                                                                                |   |      |              |
|      | Feedstuff, feed, grain 2 µg/kg                                                                                                                                                                                                    |   |      |              |
|      | Recovery rate:                                                                                                                                                                                                                    |   |      |              |
|      | Edible oil 80-120%                                                                                                                                                                                                                |   |      |              |

Feedstuff, feed, grain 70-100%

## 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                | 1 x 96 tests |
|-------------------|--------------|
| Adhesive Strip    | 4 units      |
| Antibody          | 1 x 7ml      |
| HRP-conjugate     | 1 x 7ml      |
| Micro ELISA Plate | 1 unit       |
| Sample Diluent    | 1 x 50ml     |
| Standard          | 5 x 1ml      |
| Stop Solution     | 1 x 10ml     |
| TMB Substrate     | 1 x 12ml     |
| 10X Wash Buffer   | 1 x 50ml     |

## 相关性

The aflatoxins are a group of closely related mycotoxins that are widely distributed in nature. The most important of the group is aflatoxin B1 (AFB1), which has a range of biological activities, including acute toxicity, teratogenicity, mutagenicity and carcinogenicity. In order for AFB1 to exert its effects, it must be converted to its reactive epoxide by the action of the mixed function mono-oxygenase enzyme systems (cytochrome P450-dependent) in the tissues (in particular, the liver) of the affected animal. This epoxide is highly reactive and can form derivatives with several cellular macromolecules, including DNA, RNA and protein. Cytochrome P450 enzymes may additionally catalyse the hydroxylation (to AFQ1 and AFM1) and demethylation (to AFP1) of the parent AFB1 molecule, resulting in products less toxic than AFB1. Conjugation of AFB1 to glutathione (mediated by glutathione S-transferase) and its subsequent excretion is regarded as an important detoxification pathway in animals. Aflatoxins are well recognized as a cause of liver cancer, but they have additional important toxic effects. Aflatoxin B1 is a potent hepatocarcinogenic and mutagenic mycotoxin of *Aspergillus flavus*.

**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