

MaxSignal HTS Total Aflatoxin ELISA Kit

Description

The MaxSignal® HTS Total Aflatoxin ELISA Kit is a competitive enzyme immunoassay for the quantitative analysis of Aflatoxin B₁, B₂, M₁, G₁, and G₂. This automated solution is an end-to-end workflow platform that includes simple, comprehensive multiplexed sample prep, ready-to-use reagents and standards, and a state-of-the-art automated, high-throughput detection platform. The simplified workflow saves time, hands-on requirements, reagents and lab bench footprint and enables faster, more informed decision making around grain samples. While the solution reduces cross-contamination, additional features included in the system such as the integrated bar-code scanner can provide excellent sample traceability and can be easily linked to LIMS for seamless result recording and sharing.

DS2 Automated Laboratory ELISA

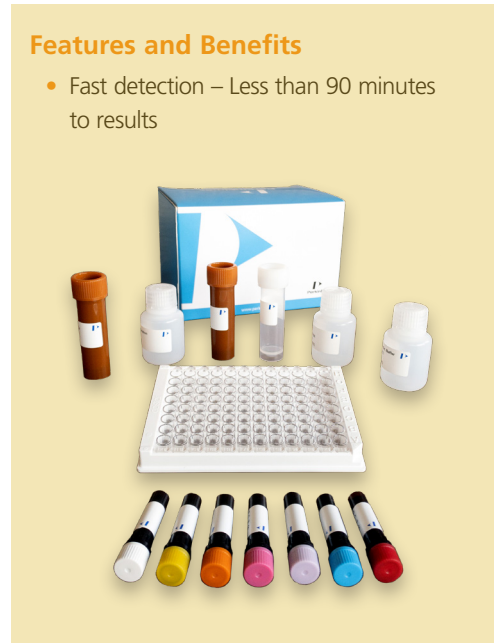
Designed with full walkaway capability, the DS2 can quickly and easily process two 96-well microplates and up to 192 samples in less than 90 minutes. Plus, it features the most intuitive control system available, including simple chain of custody and instrument diagnostics.

This system delivers no-fuss, sample-in/results-out automation of microplate assays, taking charge of the entire workflow:

- Sample dilution and distribution;
- Incubation, washing, and reagent dispensing;
- Reading with automatic data reduction and quality control;
- Automatic barcode scanning.

Features and Benefits

- Fast detection – Less than 90 minutes to results



| CATALOG # | PRODUCT | QTY |
|--------------|--|----------|
| FOOD-4008-01 | MaxSignal® HTS Total Aflatoxin ELISA Kit | 96 tests |
| DS2010 | DS2 Automated Laboratory ELISA | |

For more information visit www.perkinelmer.com

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2020 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

FOOD_SHT_20201102_1.1 PKI