

MaxSignal HTS Automated Seafood Solution

Description

Designed and developed specifically for the aquaculture industry, the PerkinElmer MaxSignal® HTS Nitrofurans and Chloramphenicol ELISA Kits and DS2 Automated Laboratory ELISA solution delivers a simple, simultaneous 5-in-1 sample preparation method for AOZ, AMOZ, SEM and AHD Nitrofurans as well as Chloramphenicol. The simplified workflow saves time, hands-on requirements, reagents and lab bench footprint and enables faster, more informed decision making around incoming seafood lots. 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

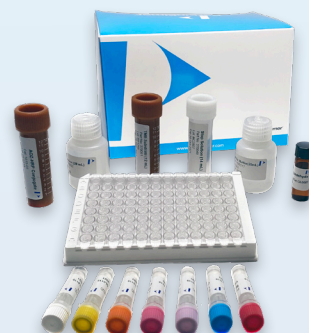
Designed with full walkaway capability, the DS2 can quickly and easily process two 96-well microplates and up to 192 samples in less than 90 minutes. Plus, it features the most intuitive control system available, including simple chain of custody and instrument diagnostics.

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

- Save money – 5-in-1 method requires one-fifth of the organic solvents for extraction
- Faster detection – Less than 90 mins from sample extract to results
- Easy, highly accurate sample preparation



| CATALOG # | PRODUCT | QTY |
|--------------|---|----------|
| FOOD-4001-01 | MaxSignal® HTS Chloramphenicol ELISA Kit | 96 tests |
| FOOD-4002-01 | MaxSignal® HTS Furazolidone (AOZ) ELISA Kit | 96 tests |
| FOOD-4003-01 | MaxSignal® HTS Furaltadone (AMOZ) ELISA Kit | 96 tests |
| FOOD-4004-01 | MaxSignal® HTS Nitrofurazone (SEM) ELISA Kit | 96 tests |
| FOOD-4005-01 | MaxSignal® HTS Nitrofurantoin (AHD) ELISA Kit | 96 tests |
| DS2010 | DS2 Automated Laboratory ELISA | |

For more information visit www.perkinelmer.com/MaxSignalHTS

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.