

Luminex[®]

ARIES[®]

Welcome to the New Way to Work



ARIES®

Welcome to the New Way to Work

The ARIES® System is crafted to increase laboratory efficiency, ensure result accuracy, and fit seamlessly into today's lean laboratory.

Designed for Improved Lab Efficiency and Workflow

- Universal Assay Protocol may enable multiple sample types and up to 12 different assays to be run together in a Random Batch.
- ARIES® software and user interface allow you to control every aspect of the sample to answer environment in your laboratory.
- Two independent modules allow for simultaneous STAT and batch testing on a single ARIES® System.
- Electronic reporting accelerates delivery of results.

Designed for Seamless Integration and Testing Accuracy

- Internal barcode scanning matches samples to cassettes, enabling Position Independent Results no matter where each cassette is placed.
- Integrated sample processing control ensures the assay run is successful from extraction through amplification.
- Bidirectional LIS connectivity enables electronic ordering.*
- Integrated touchscreen PC eliminates the need for computer peripherals, maximizing valuable bench space.

*Validation of the LIS compatibility must be performed by the user.

Workflow



Step 1
Load the sample
into the cassette



Step 2
Insert up to six cassettes
into the magazine



Step 3
Insert the magazines
into the instrument



Step 4
Review &
release results

Specifications

Physical Dimensions (Approximate)	39.1 cm x 61.2 cm x 95.0 cm (15.4" x 24.1" x 37.4")						
Weight	63.5 kg (140 lb)						
Power Supply	120 volts (US) or 220 volts (EU)						
Startup Time	Less than 5 minutes						
Hands-on Time**	Less than 2 minutes/sample						
Time to Results	Less than 2 hours						
Throughput	12 samples in less than 2 hours 48 results in 8 hours						
Sample Input Volume	200-400 µL						
Processing	2 Independent modules (1-6 samples/module); allows Random Batch and STAT testing						
Data Input	Handheld/internal bar scanning/manual						
PC	Integrated touchscreen						
Operating System	Microsoft® Windows Embedded 7						
Consumables	ARIES® assay cassettes						
Ancillary Equipment (optional)	Printer for ARIES® instrument						
Heat Source	Peltier						
Heat Output (BTU)	2730						
Operating Temperature	15° C to 35° C (59° F to 95° F)						
Temperature Control	55° C to 95° C (131° F to 203° F) within +/- 1.0° C						
Connectivity	Network and 5 x USB Type-A ports						
Network Settings	Integrates bidirectionally with LIS either through HL7 or CSV file formats						
Detector	Photodiode						
Optical Channels	Channels	1	2	3	4	5	6
	Ex/Em (nm)	390/475	464/517	525/556	557/588	594/624	630/690

**Some assays may require preparation prior to loading the cassette.

Ordering Information

Product Name	
Product Name	ARIES® Two Module System (IVD)
Part Number	
Part Number	ARIES-M12V1-IVD
ARIES® System Core Components Included	
Instrument	
Two Magazines	
Two Sample Prep Trays	
Handheld Barcode Scanner and Stand	
System Operation Manual IVD	
Quick Guide IVD	
Power Cord	
Optional Components	
Printer for ARIES® Instrument, Dell C1760nw	CN-PC13-01 (US) For all others, contact your country office for information.



orders@luminexcorp.com or support@luminexcorp.com

For In Vitro Diagnostic Use. Products are region specific and may not be approved in some countries/regions. Please contact Luminex at support@luminexcorp.com to obtain the appropriate product information for your country of residence. The ARIES® Systems are class 1(I) laser products.

©2015-2017 Luminex Corporation. All rights reserved. Luminex and ARIES are trademarks of Luminex Corporation, registered in the U.S. and other countries. Microsoft Windows is a trademark of Microsoft Corporation.

www.luminexcorp.com

HEADQUARTERS	EUROPE	CANADA	CHINA	JAPAN
UNITED STATES				
+1.512.219.8020	+31.73.800.1900	+1.416.593.4323	+86.21.802.31150	+81.3.5545.7440
info@luminexcorp.com	europe@luminexcorp.com	info@luminexcorp.com	info@luminexcorp.com	infojp@luminexcorp.com