

Luminex

ARIES[®] M1

Welcome to the New Way to Work



A New Addition to the ARIES® Sample to Answer System

The ARIES® M1 System* shares the same cassette-based sample to answer molecular diagnostic workflow as the ARIES® System, reducing hands-on time and simplifying operations. Like its predecessor, the ARIES® M1 System also shares the following:

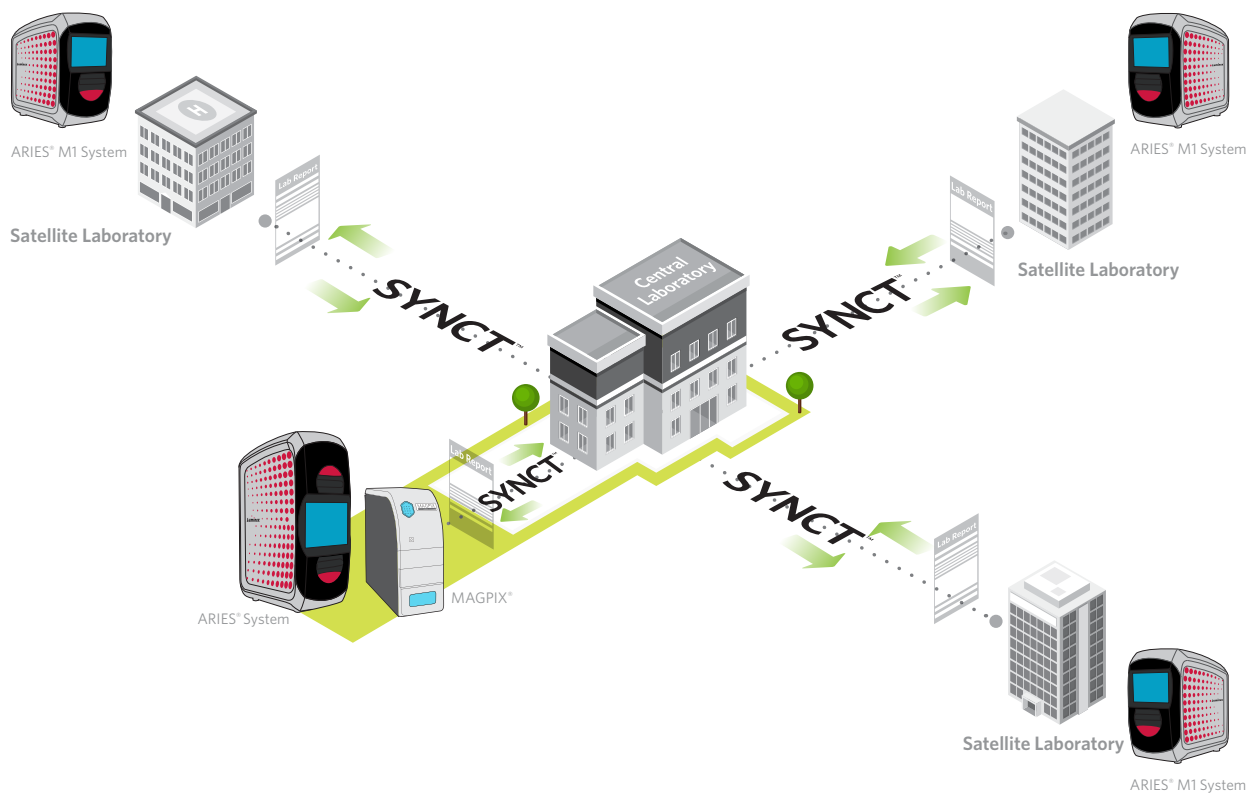
- Ability to run same IVD Assay menu as the ARIES® System
- Universal Assay Protocol may enable multiple sample types, including laboratory developed tests (LDTs), and up to six different IVD assays to be run together in a Random Batch
- Control over all aspects of the sample to answer environment through the ARIES® software and user interface
- Bidirectional laboratory information system (LIS) connectivity
- Position Independent Results based on internal barcode scanning that matches samples to cassettes
- Strong process quality from extraction through amplification with integrated sample processing controls

** Product is FDA Cleared, CE Marked, and Health Canada Licensed for IVD use.*



Enable Molecular Testing in Your Satellite Lab

The ARIES® M1 System's affordable price point enables all laboratories within a hospital network, including smaller satellite laboratories with lower throughput, to standardize molecular testing. Bidirectional laboratory information system communications and SYNCT™ Software send results instantly to the central laboratory, bringing together a suite of Luminex system solutions.



Added Flexibility in Implementing Molecular Diagnostics

Accommodate critical testing demands by supplementing your ARIES® System with an ARIES® M1 System. The ARIES® M1 System can be used for additional batched throughput or can be reserved for STAT testing requirements that require immediate processing.

Specifications

Physical Dimensions (Approximate)	39.1 cm x 61.2 cm x 64.5 cm (15.4" x 24.1" x 25.4")						
Weight	35.6 kg (78.5 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	6 samples in less than 2 hours 24 results in 8 hours						
Sample Input Volume	200-400 µL						
Processing	One module (1-6 samples/module)						
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)	850						
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.

Luminex
complexity simplified.

To learn more, please visit: www.luminexcorp.com/ARIES

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. ARIES Systems are class 1(I) laser products. Validation of the LIS compatibility must be performed by the end user.

©2016-2017 Luminex Corporation. All rights reserved. Luminex, MAGPIX, and ARIES are trademarks of Luminex Corporation, registered in the U.S. and other countries. SYNCT is a trademark of Luminex Corporation.

HEADQUARTERS

UNITED STATES

+1.512.219.8020

info@luminexcorp.com

EUROPE

+31.73.800.1900

europe@luminexcorp.com

CANADA

+1.416.593.4323

info@luminexcorp.com

CHINA

+86.21.802.31150

info@luminexcorp.com

JAPAN

+81.3.5545.7440

info@luminexcorp.com

BR19154