

## ARIES® MRSA Assay



The ARIES® MRSA Assay is a rapid, sample to answer, PCR-based *in vitro* diagnostic test for the direct detection of methicillin-resistant *Staphylococcus aureus* (MRSA) DNA from nasal swabs, and is intended as an aid in the prevention and control of MRSA infections.

- **Clinically Relevant Targets** — Incorporation of *mecA/mecC*, *SCCmec*, and *orfX* for a single MRSA call
- **Streamlined Workflow** — On-board sample processing and automated analysis make running the assay quick and easy
- **Rapid Turnaround Time** — Results are available in less than two hours
- **Flexibility and Efficiency** — Run up to 12 ARIES® assays simultaneously, including ARIES® MRSA, *C. difficile*, and Group B Strep
- **Built-In Patient Safety** — Internal barcode scanning matches samples to cassettes, decreasing the risk of sample input errors

### Workflow

The operator transfers 300 µL of sample from the ESwab™ container to the Sample Processing Tube, vortexes, and transfers 200 µL of processed specimen to the cassette. The cassette is then placed into an ARIES® Magazine, loaded into the ARIES® System, and the run is initiated.



**Step 1**  
Prepare the sample using the provided Sample Processing Tube



**Step 2**  
Load the sample into the cassette



**Step 3**  
Insert up to six cassettes into the magazine



**Step 4**  
Insert the magazines into the instrument



**Step 5**  
Review & release results

**Table 1. ARIES® MRSA Assay Performance Compared to Direct and Enriched Culture**

	Sensitivity	Specificity
<b>ARIES® MRSA Assay</b>	97/104 (93.3%)	1,550/1,658 (93.5%) <sup>1</sup>

Refer to Package Insert for additional details: Luminex Corporation | ARIES® MRSA Assay Kit Package Insert.

## Performance

Performance of the ARIES® MRSA Assay was evaluated prospectively at four geographically distinct clinical sites within the United States using clinical nasal swab samples.

## Usage and Targets

To ensure reliable identification of MRSA, the ARIES® MRSA Assay detects several targets:

- **mecA/mecC** — These genes have been shown to cause high-level resistance to methicillin by coding for alternative penicillin-binding proteins.
- **SCCmec** — SCCmec is a mobile genetic element, or cassette, that can carry *mecA/mecC*. The ARIES® MRSA Assay targets the region between the *Staphylococcus aureus* genome and this cassette to determine if the *mec* gene is specifically found in *S. aureus*.
- **orfX** — The C terminus of this gene is used to aid in identification of *S. aureus*.

The ARIES® MRSA Assay is designed to identify MRSA from nasal swab specimens collected using the Liquid Amies Elution Swab (ESwab™) Collection and Transport System, or equivalent. The ARIES® MRSA Assay is indicated for use with ARIES® Systems.

## Ordering Information

Product Name	Part Number
<b>ARIES® MRSA Assay Kit</b> Includes: 24 ARIES® MRSA Cassettes                      25 ARIES® Sample Processing Tubes	50-10045
<b>ARIES® MRSA Sample Processing Tubes (25)</b>	11-00516
<b>ARIES® MRSA Assay Protocol File Kit</b>	CN-0384-01
<b>ARIES® Two Module System</b> Includes: Instrument Two Sample Prep Trays Two Magazines System Operation Manual Quick Guide Handheld Barcode Scanner and Stand Power Cord	ARIES-M12V1-IVD
<b>ARIES® M1 System</b> Includes: Instrument Two Sample Prep Trays One Magazine System Operation Manual Quick Guide Handheld Barcode Scanner and Stand Power Cord	ARIES-M6V1-IVD
<b>SYNCT™ Software</b>	CN-SW47

Products are FDA cleared for IVD use.



**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.

©2019 Luminex Corporation. All rights reserved. Luminex and ARIES are trademarks of Luminex Corporation, registered in the U.S. and other countries. SYNCT is a trademark of Luminex Corporation. ESwab is a trademark of COPAN Diagnostics Inc.

[www.luminexcorp.com](http://www.luminexcorp.com)

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

+86.21.8036.9888

info@cn.luminexcorp.com

JAPAN

+81.3.5545.7440

info@jp.luminexcorp.com

55205554