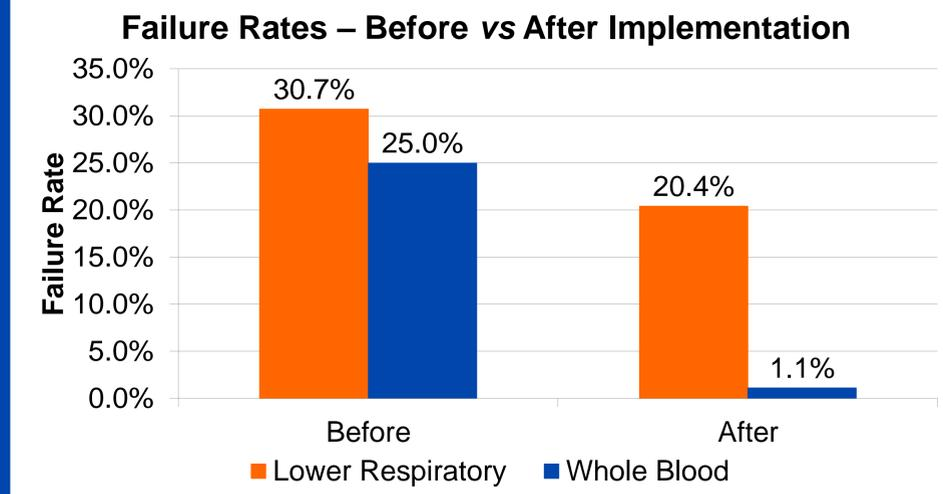
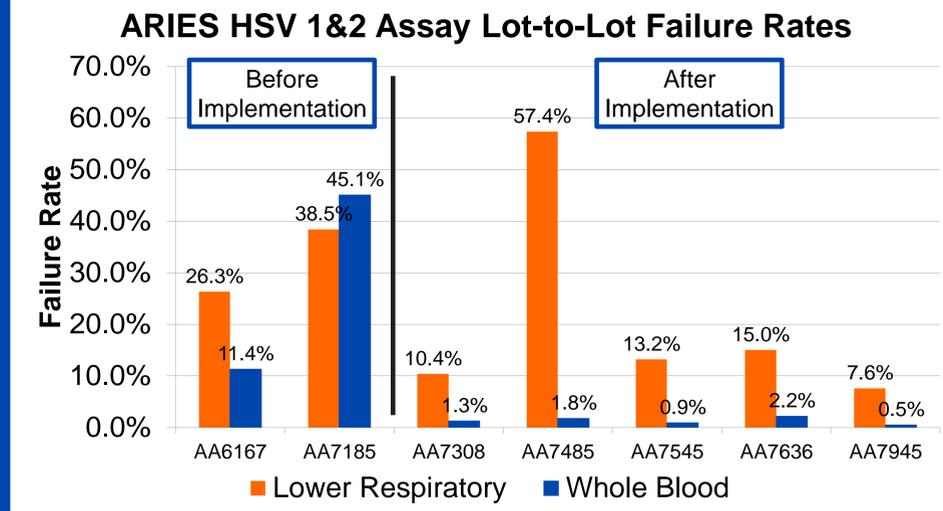


Implementation of a viscous fluid protocol led to a significant decrease in invalid rates for both whole blood and lower respiratory specimens tested on the Luminex ARIES.

Savings (time and \$\$\$) ↑
 ↓ Turnaround time

- One lower respiratory specimen yielded a negative result prior to treatment, but was positive for HSV-2 following treatment with the viscous fluid protocol. The untreated specimen was positive on repeat testing.
- One lower respiratory specimen was invalid following the viscous fluid protocol, but yielded a positive HSV-1 result on repeat testing.



CONCLUSIONS

- Pretreatment of lower respiratory specimens using a viscous fluid protocol showed similar assay performance compared to non-treated specimens, with treated whole blood specimens exhibiting increased assay sensitivity compared to non-treated specimens.
- The LoD in both specimen types remained the same for both HSV-1 and HSV-2 following the viscous fluid protocol.
- Implementation of this process led to a significant decrease in the invalid rates for both specimen types, resulting in cost-savings for the laboratory and decreased assay turnaround time.

Validation of a Viscous Fluid Preparation Protocol for Treating Whole Blood and Lower Respiratory Specimens Prior to Testing by the Luminex ARIES HSV 1&2 Assay

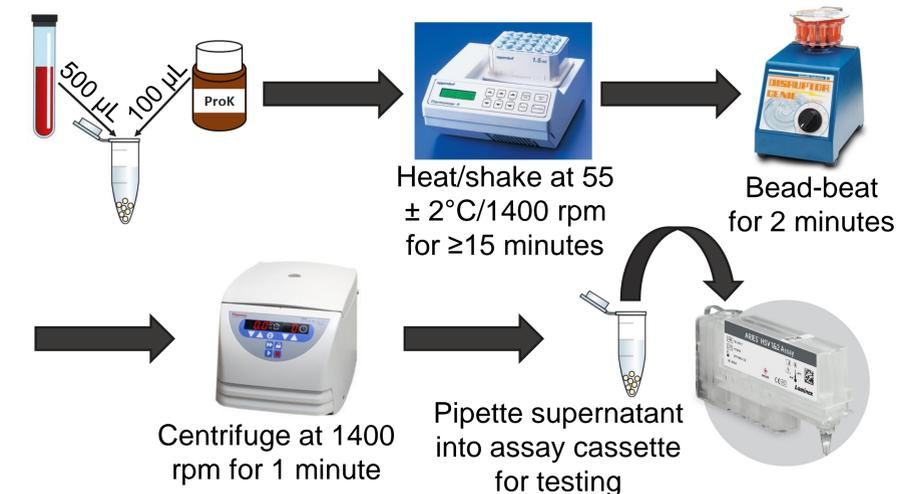
A. C. Boerger, B. J. Duresko, and M. J. Binnicker
 Mayo Clinic, Rochester, MN

BACKGROUND

- Diagnosis of herpes simplex virus (HSV) is routinely made based on clinical findings and supported by laboratory testing using polymerase chain reaction (PCR) or viral culture.
- The ARIES HSV 1&2 assay (Luminex) is a qualitative, real-time PCR test for the direct detection and differentiation of HSV-1/2 that is FDA-cleared for testing cutaneous and mucocutaneous specimens collected from symptomatic patients.

METHODS

- Following implementation of the ARIES HSV 1&2 assay, high rates of invalid results were observed for two off-label, validated sources (38.5% for lower respiratory and 45.1% for whole blood).
- Six analyte-negative lower respiratory and whole blood specimens were spiked with positive control material (ZeptoMetrix) at the previously established limit of detection (LoD) for each respective analyte and source.
- These specimens were treated using the protocol illustrated below and tested by the ARIES.



RESULTS

| Average Change (Ct Value) Before vs After Treatment | | |
|---|-------|-------|
| Specimen Source | HSV-1 | HSV-2 |
| Lower Respiratory | +0.1 | 0.0 |
| Whole Blood | -1.8 | -1.2 |



Scan the QR code to download the full abstract

