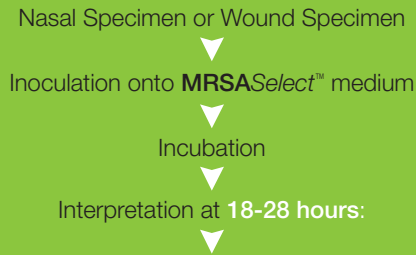


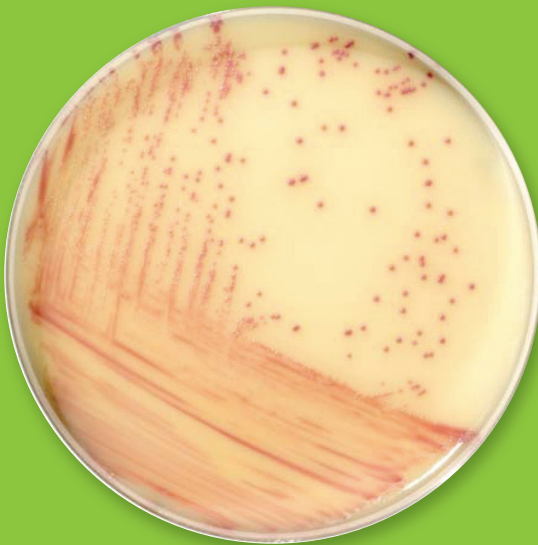
## MRSASelect™ Procedure and Interpretation



Positive

OR

Negative



*Pink colonies = MRSA*

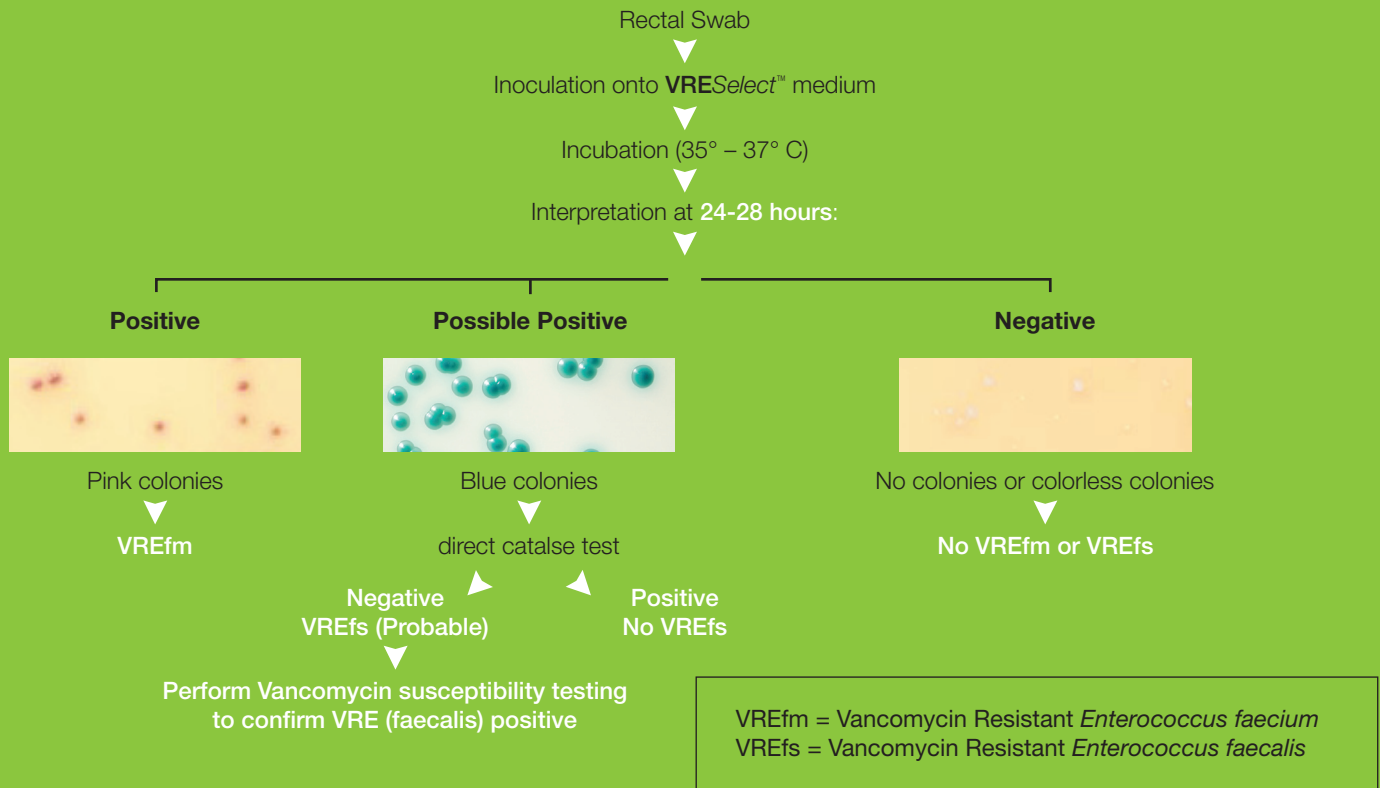


*No Growth or White colonies = No MRSA*

 VRESelect™ on reverse side

**BIO-RAD**

## VRESelect™ Procedure and Interpretation



 MRSASelect™ on reverse side

For more information, visit us on the web at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics)  
Call toll free at 1-800-2BIO-RAD (1-800-224-6723, option 2, option 3);  
Outside the US, contact your local sales office.

