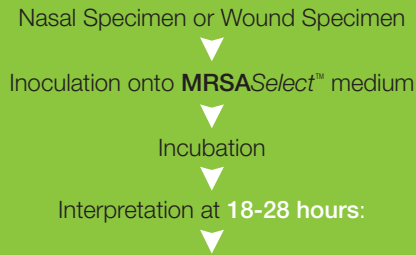


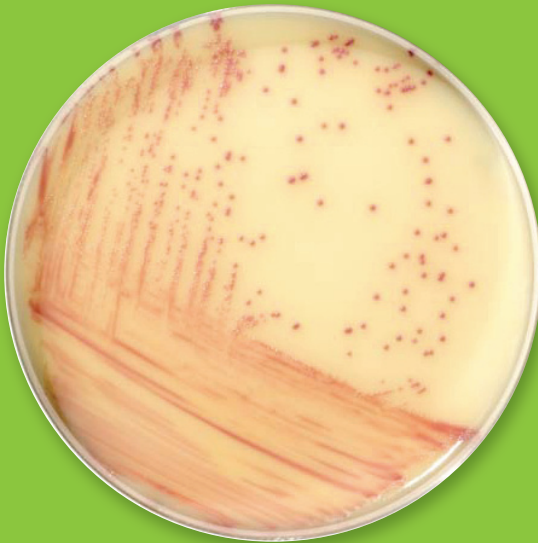
MRSASelect™ Procedure and Interpretation



Positive

OR

Negative



Pink colonies = MRSA

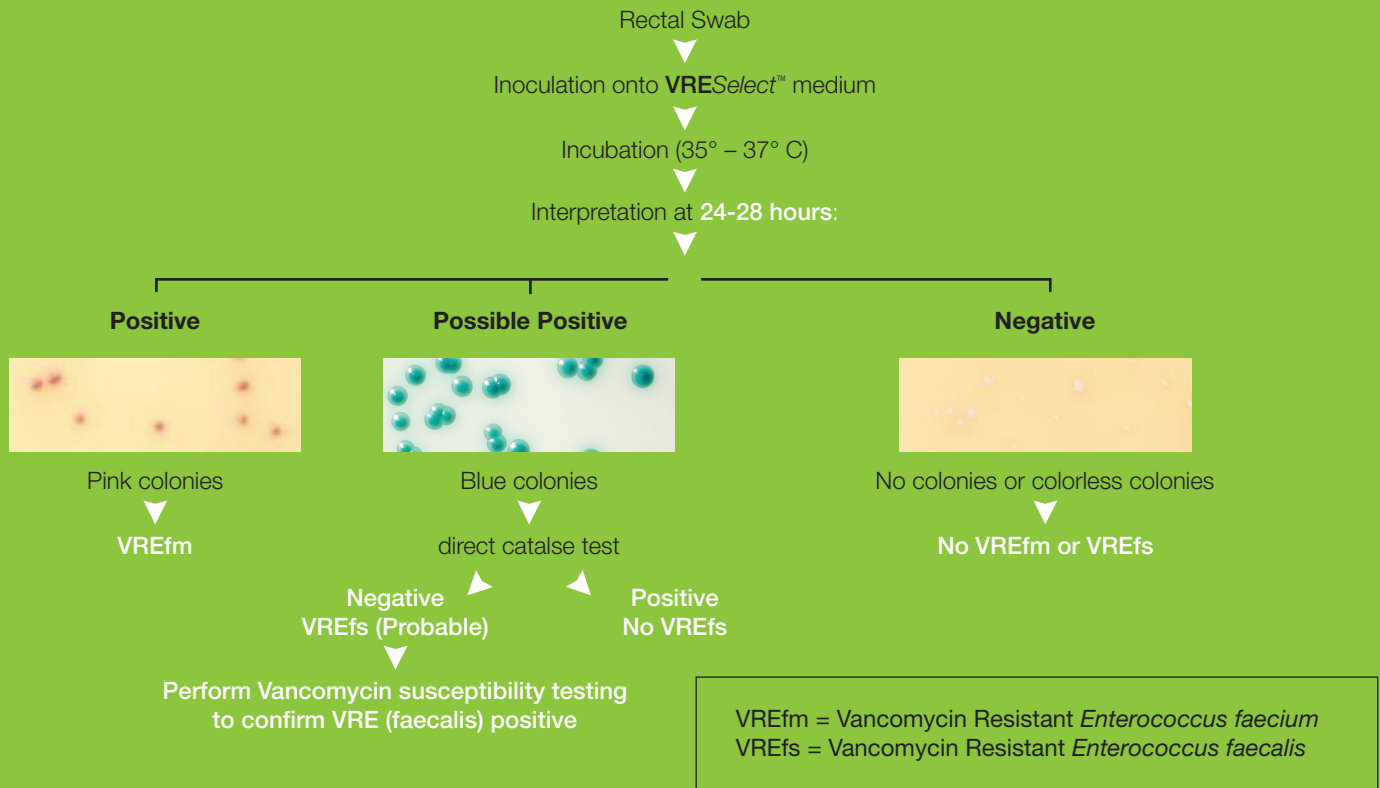


No Growth or White colonies = No MRSA

 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
Call toll free at 1-800-2BIO-RAD (1-800-224-6723, option 2, option 3);
Outside the US, contact your local sales office.

