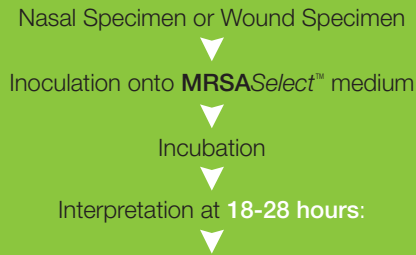


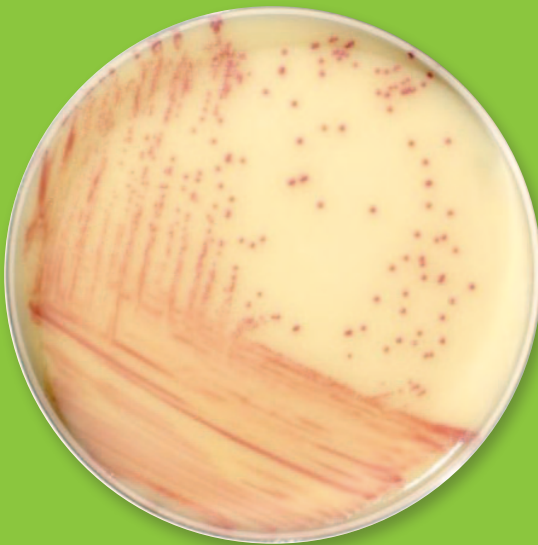
MRSASelect™ Procedure and Interpretation



Positive

OR

Negative



Pink colonies = MRSA



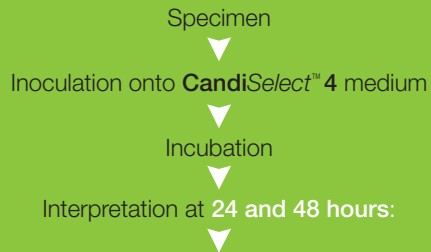
No Growth or White colonies = No MRSA



CandiSelect™ 4 on reverse side

BIO-RAD

CandiSelect™ 4 Procedure and Interpretation



Direct Identification



Candida albicans
Pink to purple colored colonies

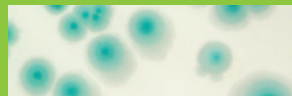
Presumptive Identification



Candida tropicalis
Intense turquoise colored colonies, spherical,
regular outline – smooth morphology



Candida glabrata
Turquoise colonies, shiny, flat,
regular outline – smooth morphology



Candida krusei
Large turquoise colonies, dry appearance,
irregular outline – rough morphology



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.

BIO-RAD