



# MRSASelect™

## The Direct Identification

### Optimal Selectivity

- Selective detection and direct identification of MRSA in **18 – 28** hours
- High sensitivity and specificity

### Ease-of-Use

- Simple procedure: inoculation, incubation and reading
- Easy to read and interpret results

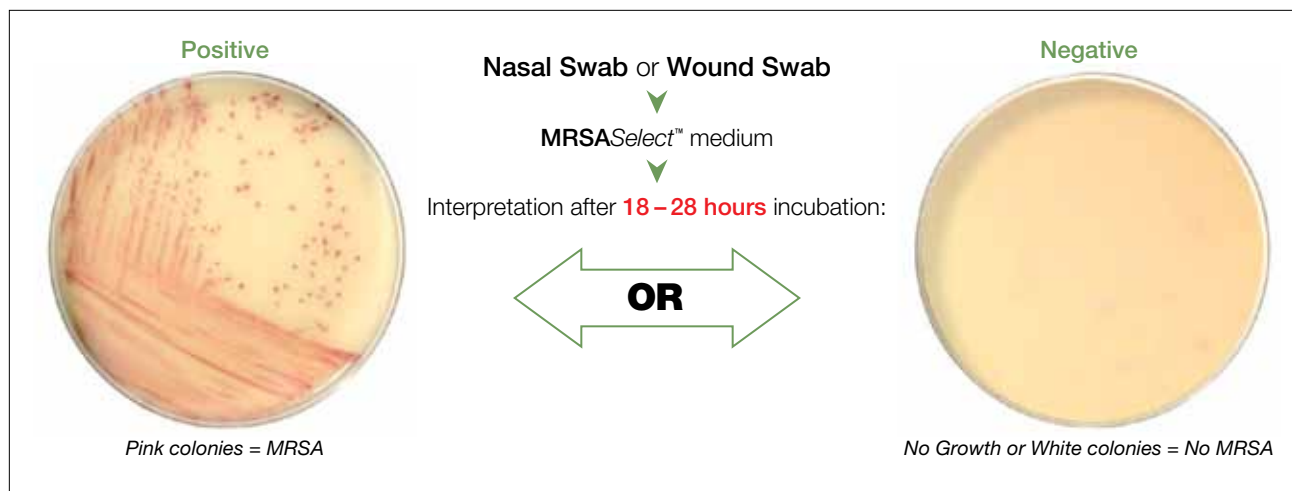
### **WOUND SPECIMEN CLAIM**

Now FDA-cleared for use with both nasal and **WOUND** specimen types



**BIO-RAD**

**MRSASelect™ Procedure**



**Performance Characteristics**

The performance of MRSASelect™ with anterior nares specimens was compared to a routine culture and to a commercially available chromogenic medium.

MRSASelect™ 24 hours	Routine Culture			
	Pos	Neg	Total	% Agreement
<b>Pos</b>	227	33	260	96%
<b>Neg</b>	10	1502	1512	98%
<b>Total</b>	237	1535	1772	

MRSASelect™ 24 hours	Commercial Chromogenic Media – 48 hours			
	Pos	Neg	Total	% Agreement
<b>Pos</b>	297	24	321	94%
<b>Neg</b>	18	2674	2692	99%
<b>Total</b>	315	2698	3013	

The performance of MRSASelect™ with skin and soft tissue wound specimens was compared to a routine culture and to TSB.

MRSASelect™	Routine Culture and TSB			
	Pos	Neg	Total	% Agreement
<b>Pos</b>	209	4	213	92%
<b>Neg</b>	19*	711	730	99%
<b>Total</b>	228	715	943	

\* For 12 out of 19 samples, MRSA was isolated only from TSB with 6.5% NaCl and were not isolated on initial direct culture.

**Note:** These tables show percent agreement and not sensitivity.

Further testing was performed to extend the incubation time

MRSASelect™ Results	Incubation Times				% Agreement
	18h	20h	24h	28h	
<b>No. Positive Samples</b>	21	21	21	21	100%
<b>No. Negative Samples</b>	179	179	179	179	100%

**Ordering Information**

**Catalog No.**      **Description**  
63747      MRSASelect™ ..... 20 plates



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics)*

Clinical  
Diagnostics Group

**Website** [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400 **Canada** 1-514-334-4372 **China** 86-21-64260808 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00 **Germany** +49-(0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 1-800-180-1224 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** +52 (55) 5488-7670 **The Netherlands** +31-318-540666 **New Zealand** 64-9-415-2280 **Norway** 47-23-38-41-30 **Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000