

---

## Spectra™ MRSA Evaluation Panel

---

Spectra™ MRSA Evaluation Panel is comprised of ready-to-use, disposable inoculating loops containing stabilized, viable microorganisms. This Laboratory Use product is formulated for use to demonstrate colony morphology and appearance on Spectra™ MRSA (REF R01821) surface. The microorganisms contained in each loop should be inoculated onto the agar surface according to the following instructions:

### Direct Streak Method

**Note:** Recommended when inoculating a single type of media.

1. Warm the Spectra™ MRSA plate to room temperature.
2. Remove the red sheath from the loop.
3. Lay the loop flat on the warm agar surface for 10-15 seconds, pressing gently onto the surface of the media. Rehydration of the film will appear as a clear "bead of moisture" upon lifting the loop from the agar surface. **Note:** The black film does not have to completely dissolve from the loop for growth of the organism to occur.
4. Streak plates for isolation and incubate according to the Spectra™ MRSA instructions for use.

### Tube Method

**Note:** Recommended when inoculating multiple types of media to avoid carryover from one type of agar to another.

1. Remove the red sheath from the loop. Break the loop shaft off from the handle directly into a tube containing 0.5-1.0 ml of liquid medium (e.g., Tryptic Soy Broth).
2. Incubate the tube at 35-37°C for 10-20 minutes to ensure the black film has completely dissolved from the loop.
3. Shake the tube gently to resuspend the organism.
4. Inoculate appropriate media with a transfer pipette. Multiple plates may be inoculated at this time.
5. Streak plates for isolation, incubate, and follow Spectra™ MRSA instructions for use.

LC4847

2011-02-15

**remel**

12076 Santa Fe Drive, Lenexa, KS 66215, USA

General Information: (800) 255-6730 Website: [www.remel.com](http://www.remel.com) Email: [remel@remel.com](mailto:remel@remel.com)

Local/International Phone: (913) 888-0939 International Fax: (913) 895-4128