

## Spectra™ VRE Evaluation Panel

Remel Spectra™ VRE 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™ VRE (REF R01830) agar plates. The microorganisms contained in each loop should be inoculated directly onto the agar plate according to the following instructions:

### Direct Streak Method

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

1. Warm the Spectra™ VRE 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™ VRE instructions for use.

### Tube Method

**Note:** Recommended when inoculating multiple types of media to avoid carryover of 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. Shake the tube gently to resuspend the organism. Inoculate appropriate media with a transfer pipette. Multiple plates may be inoculated at this time.
3. Streak plates for isolation, incubate, and follow Spectra™ VRE instructions for use.