

**VCF 250/Additive****355-5532****DEFINITION**

The VCF mixture (Vancomycin-Colimycin-Fungizone) is used as a selective inhibitor antibiotic in media used for the culture of *Neisseria*:

- Gonococci-Meningococci medium + Supplement G + VCF 250

**PRESENTATION****VCF 250**

5 ml x 1 vial

**code 355-5532****STORAGE**

- Ready to use: + 2°C to 8°C.
- Expiration date and batch number are shown on the package.

**QUALITY CONTROL OF MANUFACTURER**

Every product manufactured and marketed by Bio-Rad is subject to a quality-assurance procedure at all stages, from the reception of raw materials to the marketing of the end-product. Each batch of finished product undergoes quality control and is marketed only if it satisfies the acceptability criteria.

Documentation relative to the production and control of each batch is kept on file.