

Multivitamin Supplement

355-5592

DEFINITION

The multivitamin supplement is used to enrich media used for the culture of fastidious bacteria, *Neisseria* and *Haemophilus* in particular.

PRESENTATION

Ready to use

1 ml x 6 vial

code 355-5592

STORAGE

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

THEORETICAL FORMULA AND UTILISATION

See the technical sheet for "Chocolate agar + multivitamin supplement"

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.