



Globular Extract

355-4241

DEFINITION

Globular extract is prepared from defibrinated horse blood. It is separated by heating, then sieving and filtration.

PRESENTATION

Ready to use

10 ml x 1 vial

code 355-4241

STORAGE

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

UTILISATION

Globular extract contains growth factors X and V. When added to the usual liquid media, it permits abundant culture of bacteria with difficult growth.

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.