

## Supplement G

356-2692

### DOMAINE D'APPLICATION

Supplement G, the components of which are used as growth factors, make it possible to obtain a medium suitable for the culture of *Neisseria* and *Haemophilus*.

### PRESENTATION

#### Ready to use

50 ml x 1 bottle

code 356-2692

### STORAGE

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

### THEORETICAL FORMULA

Horse plasma	330 ml
Yeast extract	200 ml
Globular extract	470 ml
Glucose	1.3 g

### 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.