

Meat-Liver Glucose/Broth

355-3614

DEFINITION

Non-selective liquid medium conducive to the growth of numerous aerobic and anaerobic bacteria.

PRINCIPLE

The nutrient substances provided by the meat-liver base, and the glucose used as an energy source, favor the growth of most anaerobic bacteria.

PRESENTATION

Ready-to-use

10 ml x 25 tubes

code 355-3614

STORAGE

- Ready-to-use: + 2°C to 8°C.
- Dehydrated: + 15°C to 25°C, in carefully-sealed bottles in a cool, dry place.
- Expiration date and batch number are shown on the package.

THEORETICAL FORMULA

Meat-liver base	29.5 g
Glucose	2 g
Cysteine hydrochloride	0.5 g
Distilled water	1,000 ml
Final pH (25°C) = 7.4 ± 0.2	

OTHER PRODUCTS REQUIRED (NOT SUPPLIED)

- Diluent(s)
- Distilled water

EQUIPMENT REQUIRED (NOT SUPPLIED) (non-exhaustive)

- Scales
- Sterile weighing bags
- Grinder
- Hotplate
- Mixer-homogenizer
- Test tubes (16 x 160 mm) with autoclave-proof stoppers and corresponding Durham Bells
- Sterile pipettes (code 355-0751) or inoculating loop
- Water-bath precise to ± 1°C
- Thermostatically-controlled incubator or incubation room, precise to ± 1°C
- Autoclave
- All usual laboratory equipment.

PROTOCOL

Introduce 1 ml or an inoculating loop of the specimen to be analyzed into a tube of regenerated broth cooled to 44°C - 47°C.

Incubate at 37°C for 24 - 72 hours.

Then, using an inoculating loop, isolate on a suitable for the test being carried out.

PRECAUTIONS

- The time lapse between the end of the preparation of the stock solution (or the 10⁻¹ dilution in the case of a solid product) and the moment when the dilutions come into contact with the culture medium must not exceed 15 minutes.
- To avoid re-oxygenation of the medium, care should be taken not to shake it violently.
- Comply with Good Laboratory Practice.

PERFORMANCES / QUALITY CONTROL OF THE TEST

The growth performances of the media are verified with the following strains:

STRAINS	Results after 24 - 72h culture at 37°C
<i>Clostridium perfringens</i> ATCC 13124	Good growth
<i>Peptostreptococcus anaerobius</i> CIP 602	Good growth
<i>Clostridium sporogenes</i> ATCC 19404	Good 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.

KEY WORDS

Meat-liver-Glucose / Aerobic and anaerobic bacteria / Growth / Glucose / Non-selective / Medium.