

Sabouraud/Broth

355-6614
356-4854

DEFINITION

Sabouraud liquid medium is recommended by the French Pharmacopeia Code for carrying out sterility tests on pharmaceutical products. This medium permits the culture of yeasts, molds and acidophilic bacteria.

PRINCIPLE

The presence of the nutrient substances provided by meat peptic peptone and tryptic casein hydrolysate favors the growth of yeasts and molds. The glucose is used as an energy source.

PRESENTATION

• Ready-to-use

10 ml x 25 tubes

code 355-6614

• Dehydrated

500 g

code 356-4854

THEORETICAL FORMULA

| | |
|-----------------|-----------|
| Peptone | 11 g |
| Glucose | 20 g |
| Distilled water | 1,000 ml |
| Final pH (25°C) | = 6.0-6.3 |

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.

OTHER PRODUCTS REQUIRED (NOT SUPPLIED)

- Distilled water

EQUIPMENT REQUIRED (NOT SUPPLIED) (non-exhaustive)

- Scales
- Sterile weighing bags
- Grinder
- Hotplate
- Mixer-homogenizer
- Test tubes (16 x 160 mm) with autoclaveproof stoppers
- 125 ml Pyrex bottles with autoclave-proof stoppers
- Sterile Pasteur pipettes (code 355-0751) or inoculating loop

- Thermostatically-controlled incubator or incubation room, precise to $\pm 1^\circ\text{C}$
- Autoclave
- All usual laboratory equipment

PREPARATION OF DEHYDRATED MEDIUM

Always shake well before use

Dissolve 30 g of powder in 1 liter of distilled water. Mix thoroughly and dispense 10 ml per tube. Sterilize in autoclave at 121°C for 20 minutes.

Reconstitution ratio: 30 g/l.

500 g of powder makes 16.6 liters of medium.

PROTOCOL

Inoculation and incubation

Inoculate the recommended quantity of the sample to be analyzed. Incubate at 20°C-25°C for 14 days (direct inoculation of medium).

READING AND INTERPRETATION

Positive tubes may present homogenous turbidity (e.g. yeasts) or remain clear with growth of molds.

PRECAUTIONS

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 of 3-10 days culture at 25°C |
|---------------------------------------|--------------------------------------|
| <i>Candida albicans</i> ATCC 26790 | Good growth |
| <i>Candida albicans</i> ATCC 10231 | Good growth |
| <i>Candida glabrata</i> | Good growth |
| <i>Candida tropicalis</i> ATCC 750 | Good growth |

Sabouraud/Broth

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

Sabouraud / Yeasts / Molds / Pharmaceutical products / Glucose / Broth / Medium.