

Agar medium C

(Sabouraud agar with glucose)
Sabouraud agar (XIX)

355-6524
355-6519
356-4494

DEFINITION

Sabouraud agar is used for the isolation, identification and culture of yeasts and molds in samples only slightly contaminated by bacteria. This medium is recommended by the French Pharmacopeia Codex for carrying out sterility test on pharmaceutical products.

Equivalent USP 30/NF 25: Medium XIX

STANDARDS

• **European Pharmacopeia 6.0** - Biological methods - 2.6.13.: Microbiological test of non-sterile products (Detection of specified micro-organisms)

• **USP 30/NF 25 US Pharmacopeia and National Formulary (2007):** Microbial Limit Tests (61) - Microbiological Tests

PRINCIPLE

The presence of the nutrient substances provided by the peptone and of the glucose, used as an energy source, favors the rapid growth of yeasts and molds. Various antibiotics can be added to make the medium inhibit bacteria.

PRESENTATION

- **Ready to use**
8 ml x 25 tubes (inclined) **code 355-6524**
100 ml x 6 bottles **code 355-6519**
- **Dehydrated**
500 g **code 356-4494**

STORAGE

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

TYPICAL FORMULA

| | |
|-----------------------------|----------|
| Peptone | 10 g |
| Glucose | 40 g |
| Agar | 15 g |
| Distilled water | 1,000 ml |
| Final pH (25°C) = 5.6 – 6.0 | |

NB: the formula has been adapted to attain the required performance criteria.

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
- 125 ml Pyrex bottles with autoclave-proof stoppers
- Sterile Petri dishes (Ø = 90 mm)
- Sterile pipettes (1 ml, etc)
- Water-bath precise to ±1°C
- Thermostatically-controlled incubator or incubation room, precise to ±1°C
- Autoclave
- All usual laboratory equipment

PREPARATION OF DEHYDRATED MEDIUM Always shake well before use.

Dissolve 42 g of powder in 1 liter of distilled water. Heat gently, shaking frequently, then bring to the boil until completely dissolved. Dispense 8 ml per tube or 100 ml per bottle and sterilize in autoclave at 121°C ± 1°C for 20 minutes.

Reconstitution ratio: 42 g/l
500 g of powder makes 11.9 liters of medium.

PROTOCOL

Inoculation and incubation

Gentamicin or Chloramphenicol can be added. Homogenize and pour into sterile Petri dishes. Using a sterile spreader, transfer the sample to be analyzed.

Incubate at 32°C ± 1°C for 3 to 7 days.

READING AND INTERPRETATION

Separately enumerate the yeasts and molds in dishes containing between 15 and 150 colonies.

Agar medium C

(Sabouraud agar with glucose
Sabouraud agar (XIX))

V3 – 12/05/11

PRECAUTION

Comply with Good Laboratory Practice.

QUALITY CONTROL

In view of the current harmonization of pharmacopeias, we recommend that you refer to the certificates of analysis for procedures relating to the quality control (performance and selectivity) of media produced by Bio-Rad.

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/Isolation/
Identification/Culture/Agar/Glucose/Medium