

Sabouraud/Agar

355-6524
355-6519
356-4494

DEFINITION

Sabouraud agar is used to isolate, identify and culture yeasts and molds in samples only slightly contaminated by bacteria. This medium is recommended by the French Pharmacopeia Code for carrying out sterility test on pharmaceutical products.

PRINCIPLE

The nutrient substances provided by the peptone, and the glucose which serves as an energy source, favor 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)
100 ml x 6 bottles

code 355-6524
code 355-6519

• Dehydrated

500 g

code 356-4494

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

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

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 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 swirling 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

Enumerate the yeasts and molds separately in dishes containing between 15 and 150 colonies.

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	Performance 24-48H at 30°C-35°C
<i>Candida albicans</i> ATCC 26790	Good growth, white
<i>Candida tropicalis</i> ATCC 750	Good growth, white
<i>Candida glabrata</i>	Good growth, white

See the second table next page...

Sabouraud/Agar

STRAINS	7 days at 30°C-35°C and 7 days at 20°C-25°C
<i>Trichophyton rubrum</i>	Good growth, downy top, back red-brown
<i>Trichophyton violaceum</i>	Good growth Violet pigment
<i>Epidermophyton floccosum</i>	Good growth, powdery, brown back light brown
<i>Microsporum canis</i>	Good growth, downy, back yellow-orange

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