

Sabouraud + Chloramphenicol/Agar

355-6599
356-4644

DEFINITION

Sabouraud Chloramphenicol Agar is recommended for the selective isolation of yeasts and filamentous fungi (Dermatophytes and other fungi) from biological specimens presenting mixed fungal and bacterial flora.

PRESENTATION

Ready to use

100 ml x 6 bottles

code 355-6599

• Dehydrated

500 g

code 356-4644

THEORETICAL FORMULA

Peptone	10 g
Glucose	20 g
Chloramphenicol	0.2 g
Agar	15 g
Distilled water	1,000 ml
Final pH (25°C) = 6.0 ± 0.2	

STORAGE

- Ready to use: +2-25 °C
- Expiration date and batch number are shown on the package.

PROTOCOL

Inoculation and incubation

Proceed to isolate of the specimen to be analyzed or its decimal dilutions on the Sabouraud + Actidione + Chloramphenicol agar. Incubate at 32°C for 3-7 days

PRECAUTIONS

Comply with Good Laboratory Practice.

UTILISATION

Chloramphenicol inhibits most bacterial contaminants.

PERFORMANCES/QUALITY CONTROL OF THE TEST

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

STRAINS	Performance 24-48 hr at 30-35°C
<i>Candida albicans</i> ATCC 26790	Good growth
<i>Candida tropicalis</i> ATCC 750	Good growth
<i>Candida glabrata</i>	IGood growth

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

STRAINS	24-48 hr at 35-38°C
<i>Staphylococcus aureus</i> ATCC 25923	Inhibited
<i>Pseudomonas aeruginosa</i> ATCC 27853	Inhibited

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.

Sabouraud + Chloramphenicol/Agar

Documentation relative to the production and control of each batch is kept on file.

KEY WORDS

Sabouraud + chloramphenicol / yeasts / filamentous fungi / Contaminated specimens / Medium