

## 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