

## Sabouraud + Chloramphenicol + Gentamicin/Agar

356-3774  
355-6594

### DEFINITION

Sabouraud agar with the addition of Chloramphenicol and Gentamicin is recommended for the isolation of all species of yeast and other fungi from contaminated specimens.

### PRINCIPLE

The presence of nutrient substances provided by peptone and glucose favor the growth of yeasts and moulds. Gentamicin and Chloramphenicol inhibit most bacteria.

### PRESENTATION

#### • Pre-poured

90 mm x 20 plates

code 356-3774

#### • Ready-to-use

8 ml x 25 tubes (slanted)

code 355-6594

### THEORETICAL FORMULA

Peptone	10 g
Glucose	20 g
Chloramphenicol	0.5 g
Gentamicin	0.04 g
Agar	12 g
Distilled water	1,000 ml
Final pH (25°C) = 5.8 ± 0.2	

### STORAGE

- Pre-poured: + 2°C to 20°C
- Ready-to-use: + 2°C to 25°C
- Expiration date and batch number are shown on the package.

### EQUIPMENT REQUIRED (NOT SUPPLIED) (non-exhaustive)

- Sterile pipettes (0.1 ml, etc)
- Sterile Pasteur pipettes or inoculating loop
- Sterile spreaders
- Thermostatically-controlled incubator or incubation room, precise to ± 1°C
- All usual laboratory equipment.

### PROTOCOL

#### Inoculation and incubation

Transfer 0.1 ml of sample to be analyzed or its decimal dilutions to a dish of Sabouraud + Chloramphenicol + Gentamicine medium. Spread and leave to dry. Incubate at 32°C for 3-7 days.

### READING AND INTERPRETATION

Enumerate the yeasts and molds separately and identify them by means of conventional tests and the Auxacolor® scatter.

### 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 after 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

STRAINS	Performance after 24-48H at 35°C-38°C
<i>Staphylococcus aureus</i> ATCC 25923	Inhibition
<i>Pseudomonas aeruginosa</i> ATCC 27853	Inhibition

### 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 + Chloramphenicol + Gentamicin / Yeasts / Moulds / Isolation / Glucose / Medium.