

# Sabouraud + Chloramphenicol + Gentamicin + Tetrazolium/Agar

356-3734

## DEFINITION

Medium used for the isolation of, and as an aid to, identification of yeasts from all specimens.

## PRINCIPLE

Sabouraud medium is rich and favors the rapid growth of yeasts. The selective mixture of Gentamicine and Chloramphenicol avoids bacterial contaminations liable to interfere with the reading of colonies of yeasts.

The more or less marked reduction of triphenyltetrazolium chloride enables the differentiation of certain species of yeast.

## PRESENTATION

### Ready-to-use

20 plates x 90 mm

code 356-3734

## THEORETICAL FORMULA

Peptone	10 g
Glucose	20 g
Cloramphenicol	0.5 g
Gentamicin	0.04 g
Tetrazolium	0.1 g
Agar	12 g
Distilled water	1,000 ml

Final pH (25°C) = 4.5 ± 0.2

## STORAGE

- Ready-to-use: +2°C to 8°C
- Expiration date and batch number are shown on the package.

## PROTOCOL

### Inoculation and incubation

The medium is inoculated directly with the sample. Isolation is by means of streaks using a sterile Pasteur pipette.

Incubate at 32°C for 3-7 days.

## READING AND INTERPRETATION

As soon as the colonies are well developed (24-48 hours), note the degree of Tetrazolium reduction.

## 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	Results after 24 - 48h at 30-35°C
<i>Candida albicans</i> ATCC 26790	Good growth Pinkish white colonies
<i>Candida tropicalis</i> ATCC 750	Good growth Violet colonies
<i>Candida krusei</i>	Good growth Pinkish white colonies
<i>Candida pseudotropicalis</i> CIP 513.56	Good growth Pink colonies
<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 + Gentamicine + Tetrazolium / Yeasts / Isolation / Identification / All specimens / Medium.