

Sabouraud + Actidione Chloramphenicol/Agar

355-6559

DEFINITION

Sabouraud agar with the addition of actidione and chloramphenicol is recommended for the isolation of Dermatophytes and other pathogenic fungi from heavily-contaminated specimens.

PRESENTATION

- **Ready-to-use**
100 ml x 6 bottles

code 355-6559

THEORETICAL FORMULA

Peptone	10 g
Glucose	20 g
Actidione (cycloheximide)	0.5 g
Chloramphenicol	0.5 g
Agar	15 g
Distilled water	1,000 ml
Final pH (25°C) = 6.0 ± 0.2	

STORAGE

- Ready-to-use: + 2°C to 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

Actidione inhibits the development of certain fungi (*Candida krusei*, *Candida tropicalis*, *Fusarium*, *aspergillus fumigatus*, etc) and has no action on the following pathogenic fungi, which can be isolated on this medium:

All Dermatophytes;
All *Candida* (except for *C. tropicalis*, *C. krusei*, *C. pelliculosa*);
Sporotrichum schenckii;
Histoplasma capsulatum;
Blastomyces dermatitis;
Coccidioides immitis;
Geotrichum candidum;

Trichophyton mentagrophytes;
Trichophyton rubrum;
Epidermophyton floccosum etc.

However, this medium does not allow the isolation of *Cryptococcus neoformans* or *Monosporium apiospermum*.

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

STRAINS	7 days at 30-35°C and 7 days at 20-25°C
<i>Trichophyton rubrum</i>	Good growth downy, 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
<i>Fusarium</i>	Inhibited

STRAINS	24-48H 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.

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

KEY WORDS

Sabouraud + actidione + chloramphenicol
Dermatophytes / Contaminated specimens /
Medium.