

Casitone/Agar

355-5813

DEFINITION

Casitone medium is used for examining the activity of antifungal agents using the disk-diffusion method.

Fluorocytosine must be tested on the semi-synthetic medium for fungal antibiogram.

PRESENTATION

• Ready-to-use

15 ml x 25 tubes

code 355-5813

STORAGE

• + 2°C to 8°C.

• Expiration date and batch number are shown on the package.

THEORETICAL FORMULA

Casitone	9 g
Yeast extract	5 g
Trisodium sodium citrate	10 g
Disodium phosphate	1 g
Monosodium phosphate	1 g
Charcoal-purified dextrose	10 g
Agar	20 g
Distilled water	1,000 ml
Final pH (25°C) = 6.6 ± 0.2	

PROTOCOL

The volume to be dispensed is 15 ml for a dish of 90 mm diameter.

Method: Refer to "BK Antibiogram" Technical Sheet (codes 357-0302, 357-0301, 357-0472, 357-0652).

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	Inhibition diameter (mm) after 24-48 h culture at 32°C
<i>Candida albicans</i> SDP 3153A	Nystatine: 16-25 Ampho B: 17-22 Econazole: 23-33 Clotrimazole: 18-28 Miconazole: 18-25

STRAINS	Inhibition diameter (mm) after 24-48 h culture at 32°C
<i>Candida albicans</i> SDP 41 R 5FC	Nystatine: 17-26 Ampho B: 17-22 Econazole: 25-35 Clotrimazole: 18-30 Miconazole: 20-30

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

Casitone / Disks / Antibiogram / Fluorocytosine / Medium.

BIBLIOGRAPHY

- **REGLI P., FERRARI H., GOUDARD M. et BUFFARD Y. (1982):** Intérêt du milieu Casitone pour l'étude de la sensibilité in vitro des champignons levuriformes aux antifongiques dérivés de l'imidazole, Bull. Soc. Franç. Mycol. Méd., 11: 359-362.
- **DROUHET E., et DUPOND B. (1978):** Antibiogramme des champignons aux antifongiques. Bull. Soc. Franç. Mycol. Méd.: 165-170.