

Semi-Synthetic/Agar

355-5812

DEFINITION

Semi-synthetic medium is used to study 5-fluorocytosine activity by the disk-diffusion method (in agar).

PRESENTATION

Ready-to-use

15 ml x 25 tubes

code 355-5812

STORAGE

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

THEORETICAL FORMULA

Asparagine	1.5 g
L-Histidine	10 mg
DL Methionine	20 mg
DL Tryptophan	20 mg
Ammonium sulfate	3.5 g
Purified agar	10 g
Potassium phosphate	1 g
Magnesium sulfate	0.5 g
Sodium chloride	0.1 g
Vitamins: Biotin	10 ⁻⁸ g
Thiamin	10 ⁻⁶ g
Calcium pantothenate	10 ⁻⁶ g
Pyridoxin	10 ⁻⁶ g
Nicotinic acid	10 ⁻⁶ g
Inositol	10 ⁻⁵ g
Solution of trace elements (Berthelot)	10 drops
Agar	20 g
Distilled water	1,000 ml

Final pH (25°C) = 5.6 ± 0.2

UTILISATION

The volume to be distributed is 15 ml for a plate of 90 mm diameter.

Method:

Refer to the "Antibiogram" sheet.

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

Semi-synthetic / Diffusion / Disks / Antibiogram / 5-fluorocytosine / Medium.

BIBLIOGRAPHY

DROUHET E., DUPONT B. (1978): Antibiogramme des champignons aux antifongiques, Bull. Soc. Franç. Mycol. Méd, 7, 165-170.