

Mueller-Hinton Salt Agar (+ 5 % NaCl)

355-6136

DEFINITION

Medium used to study of the susceptibility of resistant strains of *staphylococci* to oxacilline.

PRESENTATION

Ready-to-use

100 ml x 6 bottles

code 355-6136

STORAGE

- Ready-to-use: + 2°C to 8°C.

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

THEORETICAL FORMULA

Dehydrated beef infusion	4 g
Casein hydrolysate	17.5 g
Corn starch	1.5 g
Agar + 5% NaCl	10 g
Distilled water	1,000 ml
Final pH (25°C) = 7.3 ± 0.1	

EQUIPMENT REQUIRED (NOT SUPPLIED)

(non-exhaustive)

- Sterile Petri dishes (Ø = 90 mm)
- Water-bath at 100°C for fusion of ready-to-use media
- Thermostatically-controlled incubator or incubating room, precise to ± 1°C
- Oxacilline plates (code 356-6888)
- All usual laboratory equipment.

PROTOCOL

• Inoculation and incubation

Pour the melted medium, cooled to 44°C-47°C, onto Petri dishes and leave to dry for 30 minutes at 37°C.

Surface-inoculate the strain to be tested and deposit the plates of selected antibiotics.

READING AND INTERPRETATION

Reading will involve reference to our "Antibiogram" technical sheet.

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 at 37°C
<i>Staphylococcus aureus</i> CIP 107399	Resistance

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

Mueller-Hinton salt agar / *Staphylococci* / Susceptibility / Antibiogram / Oxacilline / NaCl / Medium.