

Conservation of Bacterial Strains/Agar 356-3683

DEFINITION

Medium intended mainly for the conservation of *Enterobacteria*. It can also be used for the conservation of:

- *Vibrionaceae*,
- *Pseudomonadaceae*,
- *Alcaligenes*,
- *Flavobacterium* and related bacteria,
- and possibly *Acinetobacter* and *staphylococci*.

PRINCIPLE

The nutrient substances provided by meat extract and peptone favor the conservation of bacterial strains.

PRESENTATION

Ready-to-use

50 tubes x 1 set

code 356-3683

STORAGE

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

THEORETICAL FORMULA

Meat extract	5 g
Peptone	10 g
NaCl	5 g
Agar	10 g
Distilled water	1,000 ml
Final pH (25°C) = 7.3 ± 0.2	

PROTOCOL

Inoculation is based on a culture on solid media: Nutrient Agar or Trypto-Casein-Soy Agar.

- For facultative aerobic-anaerobic bacteria: several central inoculations using a generously-loaded platinum loop.
- For aerobic bacteria, surface inoculation. Incubate at 37°C maximum for 24 hours. A bacterial strain inoculated in this medium can be kept:
 - 1 year at room temperature
 - 3 years at 4°C

Conservation is optimized if glycerol or vaseline is deposited on the surface.

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 24h culture at 37°C
<i>Staphylococcus aureus</i> ATCC 25923	Good growth
<i>Escherichia coli</i> ATCC 25922	Good growth
<i>Pseudomonas aeruginosa</i> ATCC 27853	Good growth
<i>Staphylococcus epidermidis</i> ATCC 14990	Good growth

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

Conservation of bacterial strains / *Enterobacteria* / Nutrient substances / Medium.