

## 1.3% Nutrient/Broth (with NaCl)

355-3446  
356-4065

### DEFINITION

Medium permitting the culture of non-fastidious bacteria.

### PRINCIPLE

The nutrient substances provided by peptone, meat extract and vitamin factors of yeast extract favor the growth of most bacteria.

### PRESENTATION

- **Ready to use**  
9 ml x 25 tubes **code 355-3446**
- **Dehydrated**  
500 g **code 356-4065**

### STORAGE

- Pre-poured: +2-8°C
- Dehydrated: +15-25°C, in carefully sealed bottles in a cool, dry place
- Expiration date and batch number are shown on the package.

### THEORETICAL FORMULA

Peptone	5 g
Meat extract	1 g
Yeast extract	2 g
Sodium chloride	5 g
Distilled water	1,000 ml
Final pH (25°C) = 7.4 ± 0.2	

### OTHER PRODUCTS REQUIRED (NOT SUPPLIED)

- Distilled water

### EQUIPMENT REQUIRED (NOT SUPPLIED) (non-exhaustive)

- Scales
- Sterile weighing bags
- Grinder
- Hotplate
- Mixer-homogenizer
- Vortex-type shaker
- Test tubes (16 x 160 mm) with autoclave proof stoppers
- Sterile pipettes (1 ml, etc)
- Sterile Pasteur pipettes (code 355-0751) or inoculating loop
- Thermostatically-controlled incubator or incubation room, precise to ±1°C
- Autoclave
- All usual laboratory equipment

### PREPARATION OF DEHYDRATED MEDIUM

#### Always shake well before use.

Dissolve 13 g of powder in 1 liter of distilled water. Mix well and disperse 10 ml per tube. Sterilize in autoclave at 121°C ± 1°C for 15 minutes.

**Reconstitution ratio: 13 g/l**

**500 g of powder makes 38.4 liters of medium.**

### PROTOCOL

Depending on the objective, this medium can be inoculated by:

- spreading 0.1 ml of sample on the surface,
- isolation by means of an inoculating loop or a Pasteur pipette (previously flame-sterilized).

### PRECAUTION

Comply with Good Laboratory Practice.

### PERFORMANCES/QUALITY CONTROL OF THE TEST

The growth performances of the media are verified with the following strains:

STRAINS	Result after 24 hr culture at 37°C
<i>Staphylococcus aureus</i> ATCC 25923	Good growth
<i>Staph. epidermidis</i> ATCC 12228	Good growth
<i>Salmonella Enteritidis</i> ATCC 13076	Good growth
<i>Streptococcus pyogenes</i> ATCC 19615	Good growth
<i>Escherichia coli</i> ATCC 25922	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

Nutrient1.3%/Non-fastidious/bacteria/NaCl  
/Growth/Broth/Medium