

Agar medium H with crystal violet

(MacConkey agar with crystal violet)
MacConkey agar medium with crystal violet (XVII)

356-3617
356-9084

DEFINITION

Medium used to isolate and enumerate coliform bacilli, and particularly *Escherichia coli* in the testing of microbial contamination of non-sterile Pharmacopeia products (2.6.13.).

Equivalent USP 30/NF 25: Medium XVII

STANDARDS

• **European Pharmacopeia 6.0** – Biological methods - **2.6.13.**: Microbiological test of non-sterile products (Detection of specified micro-organisms)

• **USP 30/NF 25 US Pharmacopeia and National Formulary (2007)**: Microbial Limit Tests (61) - Microbiological Tests

PRINCIPLE

The principle of the medium rests on the ability of coliforms to ferment lactose (brick-red colonies).

Due to the presence of bile salts and crystal violet, the medium inhibits Gram-positive bacteria.

PRESENTATION

• Ready to use

20 plates x 90 mm

code 356-3617

• Dehydrated

500 g

code 356-9084

STORAGE

- Ready to use: +2-20°C
- Dehydrated: +15-25°C, in carefully-sealed bottles in a cool, dry place
- Expiration date and batch number are shown on the package.

TYPICAL FORMULA

Bacteriological peptone	20 g
Bile salts	1.5 g
Sodium chloride	5 g
Lactose	10 g
Neutral red	30 mg
Crystal violet	1 mg
Agar	15 g
Distilled water	1,000 ml

Final pH (25°C) = 7.1 ± 0.2

NB: the formula has been adapted to attain the required performance criteria.

OTHER PRODUCTS REQUIRED (NOT SUPPLIED)

- Diluent(s)
- Distilled water

EQUIPMENT REQUIRED (NOT SUPPLIED) (non-exhaustive)

- Scales
- Sterile weighing bags
- Grinder
- Hotplate
- Mixer-homogenizer
- 125 ml Pyrex bottles with autoclave-proof stoppers
- Sterile Petri dishes (Ø = 90 mm)
- Sterile pipettes (1 ml, etc)
- Water-bath precise to ±1°C
- 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 51.5 g of powder in 1 liter of distilled water.

Bring to the boil until completely dissolved.

Dispense 100 ml per bottle and sterilize in autoclave at 121°C ± 1°C for 15 minutes.

Cool to 44-47°C then pour into Petri dishes.

After solidification, leave to dry in incubator at 37°C ± 1°C (partially closed).

Reconstitution ratio: 51.5 g/l

500 g of powder makes 9.7 liters of medium.

PROTOCOL

Inoculation and incubation

After presumptive screening, carry out Sub-culture on Mac Conkey medium.

Incubate at 44°C ± 1°C for 18-24 hours.

READING AND INTERPRETATION

- Lactose-negative colonies: colorless
- Lactose-positive colonies: brick-red, sometimes surrounded by an opaque halo of precipitated bile salts

Agar medium H with crystal violet

(MacConkey agar with crystal violet)
MacConkey agar medium with crystal violet (XVII)

V3 – 06/06/11

PRECAUTIONS

- The time lapse between the end of preparation of the stock solution (or the 10^{-1} dilution in the case of a solid product) and the moment when the dilutions come into contact with the culture medium must not exceed 15 minutes.
- Comply with Good Laboratory Practice.

QUALITY CONTROL

In view of the current harmonization of pharmacopeias, we recommend that you refer to the certificates of analysis for procedures relating to the quality control (performance and selectivity) of media produced by Bio-Rad.

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

MacConkey with crystal violet/Coliforms/
Escherichia coli/Food products/Isolation/
Enumeration/Lactose/Medium