

TGV Anaer/Kit

(Transport medium for live bacteria/aerobes)

355-0458

DEFINITION

TGV kit (transport of live bacteria) is a combination of swab + Amies transport medium (without charcoal). It is a soft agar.

PRINCIPLE

Amies medium is a transport medium, i.e. it ensures the survival of bacteria without promoting their multiplication.

It permits the survival of anaerobes for 24 - 48 hours at room temperature, in particular:

- *Bacteroides fragilis*,
- *Fusobacterium nucleatum*,
- *Streptococcus anaerobius*,
- *Eubacterium lentum*.

PRESENTATION

Pack of 24 kits

code 355-0458

STORAGE

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

THEORETICAL FORMULA

Sodium chloride	3 g
Potassium chloride	0.2 g
Calcium chloride	0.1 g
Magnesium chloride	0.1 g
Monopotassium phosphate	0.2 g
Disodium phosphate	1.15 g
Sodium thioglycolate	1 g
Agar	4 g
Distilled water	1,000 ml
Final pH (25°C) = 7.2 ± 0.2	

At the outset, the TGV kit consists of 2 compartments: one protecting the swab, the other containing the transport medium.

PROTOCOL

- Release the swab by removing the protective tube.
- Collect the specimen.
- Undo the swab-support lid.
- Introduce the swab into the tube containing the transport medium.
- Seal it hermetically by immersing the cotton-wool in the transport medium.

- Fill in the label:

- Product name,
- Reference,
- Collection site,
- Date,
- Time.

PERFORMANCES / QUALITY CONTROL OF THE TEST

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

STRAINS	Results at 37°C in anaerobiosis
<i>Bacteroides fragilis fragilis</i> ATCC 25285	Adequate survival over 48 hours
<i>Fusiformis fusiformis</i>	Adequate survival over 48 hours
<i>Eubacterium lentum</i>	Adequate survival over 48 hours

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

TGV / Anaerobes / Food products / Water / Transport / Medium.