

## TGV Aer/Kit

(Transport medium for live bacteria/aerobes)

355-0456

### DEFINITION

TGV (transport of live bacteria) is a combination of swab + Amies transport medium (with charcoal). It is a soft agar.

TGV can be used for the transport of strains: after collection by swab of a pure 18 - 24 hours culture, place it in the transport medium, then in special packaging for the transport of bacterial strains (**code 355-0132**).

### PRINCIPLE

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

It permits the survival of aerobes for 48 - 72 hours at room temperature, with the exception of *gonococcus* and *meningococcus*, survival of which cannot exceed 48 hours at room temperature.

Amies medium also maintains the survival of *Trichomonas vaginalis*.

### PRESENTATION

Pack of 24 sets

code 355-0456

### 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
Charcoal	10 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
<i>Neisseria meningitidis</i> ATCC 13090	Adequate survival over 48 hours
<i>Neisseria meningitidis</i> ATCC 13077	Adequate survival over 48 hours
<i>Neisseria gonorrhoeae</i>	Adequate survival over 48 hours
<i>Haemophilus influenzae</i> CIP 52151	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 / Aerobes / Food products / Water / Transport / Medium.