

MEDIUM FOR THE DIFFERENTIATION OF *ENTEROCOCCUS* SPECIES

IVD

1- INTENDED USE

Potassium tellurite agar is a medium for differentiation of *Enterococcus* species.

2- PRINCIPLE

The selectivity of this medium is due to the presence of potassium tellurite, which inhibits the growth of most bacteria associated with *Enterococcus* species.

The differentiation of *Enterococcus* species is based on their capacity to reduce potassium tellurite into tellurium, which gives the colonies a black colour.

3- HOW SUPPLIED

- Ready to use medium: Potassium Tellurite agar base
 - 25 x 20 ml tubes
- Ready to use supplement: Potassium Tellurite solution
 - 25 x 1 ml ampoules of potassium tellurite

Code 55726

4- COMPOSITION

Meat extract	8 g/l
Yeast extract	5.5 g/l
Peptone	9.5 g/l
Sodium chloride	5 g/l
Glucose	5 g/l
Agar	15 g/l

Supplement: potassium tellurite (0.8%): 1 ml/20 ml of medium.

Preparation of the medium:

Melt the agar base in a boiling water-bath, then cool to a temperature of about 50°C.

Add the contents of one 1 ml ampoule of 0.8% potassium tellurite solution to each 20 ml tube.

Mix and dispense into Petri dishes.

5- STORAGE

- Ready to use Medium/Supplement: at +2-8°C.
- The expiry date and batch number are indicated on the packaging.

6- INSTRUCTIONS**Material:**

Material provided: Potassium tellurite medium.

Inoculation:

Inoculate in very tight streaks from a fresh, pure culture on agar medium .

Incubation:

Incubate for 24 to 48 hours at 37°C.

Reading:

- Tellurite (+): black colonies (*Enterococcus faecalis*)
- Tellurite (-): no growth or white colonies (*Enterococcus faecium* or other *Enterococcus* species).

7- PERFORMANCE/QUALITY CONTROL OF THE TEST

- Appearance of the ready to use medium: **clear, amber-coloured** agar.
- Appearance of the ready to use supplement: **clear, colourless** liquid.
- The growth performances of Potassium Tellurite medium are verified with the following strains:

STRAINS	CULTURE RESULT AFTER 24-48 hours at 37°C
<i>Enterococcus faecalis</i> var <i>zymogenes</i> ATCC 29212	Black colonies
<i>Enterococcus faecalis</i> var <i>liquefaciens</i> CIP 54.30	Black colonies
<i>Enterococcus durans</i> ATCC 19432	No growth

8- QUALITY CONTROL OF THE MANUFACTURER

All manufactured reagents are prepared according to our Quality System, starting from reception of raw material to the final commercialization of the product. Each lot is submitted to quality control assessments and is only released to the market, after conforming to pre-defined acceptance criteria. The records relating to production and control of each single lot are kept within Bio-Rad.



Bio-Rad

3, boulevard Raymond Poincaré
92430 Marnes-la-Coquette France
Tel. : +33 (0) 1 47 95 60 00
Fax : +33 (0) 1 47 41 91 33



11/2003