

Sterile Distilled Water/Diluent

355-4154

DEFINITION

Sterile distilled water is generally used for the sterile reconstitution of certain products or for the standardization of inoculums (dilutions) based on samples.

STANDARDS

FOOD MICROBIOLOGY

- **NF ISO 7218/A1 (December 2001):** Food microbiology - General rules for microbiological examinations.

PRESENTATION

Ready-to-use

9 ml x 25 tubes

code 355-4154

STORAGE

- Ready-to-use: + 2°C to 25°C.
- Expiration date and batch number are shown on the package.

PERFORMANCES / QUALITY CONTROL OF THE TEST

The growth performances of the media are verified with the following strains, as a survival test:

STRAINS	Count at T0 et T45 min at 20-25°C
<i>Escherichia coli</i> ATCC 25922	± 50% col./T0
<i>Staphylococcus aureus</i> ATCC 25923	± 50% col./T0

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

Distilled water / Food products / Medium.