



Sulfamethazine Solution (0.2%) /Additive

356-2682

DEFINITION

Sulfamethazine is used in the preparation of Baird Parker medium as growth inhibitor for *Proteus*.

Its incorporation in the presentations of this medium (dehydrated and ready to use) is defined in the Baird Parker agar medium.

PRESENTATION

Ready to use

2.5 ml x 1 vial

code 356-2682

STORAGE

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

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.