

## Nalidixic Acid/Tests

**356-8618**

### DEFINITION

Nalidixic acid is used as an element of differentiation in the identification of *Campylobacter* (see "Selective supplement for *Campylobacter*")

### PREPARATION

**Disks containing 30 µg**

50 disks

**code 356-8618**

1 set of 4 cartridges

### STORAGE

- + 2°C to 8°C in a cool dry place.
- 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.