

Novobiocin/Disks

355-6350

PRESENTATION

Ready to use

30 disks at 5 µg x 1 cartridge

code 355-6350

STORAGE

- Ready-to-use: + 2°C to 8°C in a dry place.
- Expiration date and batch number are shown on the package.

PROTOCOL

Detection of susceptibility to novobiocin is carried out on Mueller-Hinton medium, following the antibiogram method (by flooding).

READING AND INTERPRETATION

Resistance to this antibiotic is defined by an inhibition diameter of less than 15 mm.

To differentiate *S. saprophyticus* from *S. cohnii* (both resistant to novobiocin), it is necessary to investigate saccharose attack under anaerobiosis: *S. saprophyticus* is positive.

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.

BIBLIOGRAPHY

- **ALMEIRA R.J., JORGENSEN J.H. (1982):** Use of Mueller-Hinton agar to determine Novobiocine susceptibility of coagulase negative *Staphylococci*, *Journal Of Clinical Microbiology*, **16**: 1155-1156.