

ONPG /Disks

355-3822

DEFINITION

Certain Enterobacteria possess:

- the enzyme necessary for the use of lactose by the bacteria (β -galactoside-permease);
- as well as the enzyme splitting the lactose molecule into glucose and galactose (β -galactosidase).

If the first of these two activities is deficient or absent, a potentially "lactose-positive" bacteria - because it possesses a β -galactosidase - would be unable to express this character and would appear "lactose-negative". Acidification of a lactose medium will manifest only later, if a reverse mutant appears, or perhaps not occur at all.

Reaction:

Orthonitrophenyl- β -D-Galactopyranoside (ONPG) in solution is colourless.

The molecule is split, like lactose, by β -galactosidase, releasing the orthonitrophenol, yellow in solution.

This reaction is carried out in buffered solution.

The disks are impregnated with ONPG + buffer.

PRESENTATION

Set of 50 disks

code 355-3822

STORAGE

- Ready to use: + 2°C to 8°C.
- Expiration date and batch number are shown on the package.
- In the event of doubts over the validity, check using a positive strain.

PROTOCOL

Bacteria, collected from agar medium, are placed in a thick (milky) suspension in 0,5 ml of distilled water in a small tube (e.g. Kahn tube). The more bacteria, the more enzymes, and therefore the reaction will be all the more rapid.

Place a disk in this suspension and incubate in a water-bath at 37°C.

READING AND INTERPRETATION

Observe after 1 hour of incubation.

The vast majority of positive reactions (yellow colour) occur in 15 to 30 minutes.

Results obtained with aerobic or anaerobic Gram-negative bacilli:

	ONPG positive release of orthonitrophenol = yellow solution	ONPG negative = colorless solution
ENTERO-BACTERIA	<p><i>E. coli</i> (the majority of strains of <i>Alkalescens dispar</i> biotype are ONPG -)</p> <p><i>Shigella sonnei</i> biotype xylose –</p> <p><i>Shigella dysenteriae</i> 1</p> <p><i>Shigella dysenteriae</i> 6</p> <p><i>Shigella boydii</i> 9</p> <p><i>Salmonella</i> sub-species III</p> <p>(<i>S. arizona</i>)</p> <p><i>Salmonella</i> sub-species II (a)</p> <p><i>Citrobacter</i></p> <p><i>Levinea malonatica</i></p> <p><i>Levinea amalonatica</i>*</p> <p><i>Klebsiella</i></p> <p>(except</p> <p><i>K. rhinoscleromatis</i>)</p> <p><i>Enterobacter Hafnia</i>*</p> <p><i>Serratia marcescens</i>*</p> <p><i>Serratia liquefaciens</i></p> <p><i>Serratia plymutica</i></p> <p><i>Yersinia</i>*</p>	<p>Other <i>Shigella</i></p> <p><i>Salmonella</i> sub-species let IV</p> <p><i>Edwardsiella</i></p> <p><i>Proteus</i></p> <p><i>Providencia</i></p>
VIBRIO-NACEAE	<p><i>Vibrio cholerae</i>*</p> <p><i>Vibrio anguillarum</i>*</p> <p><i>Plesiomonas</i></p>	<p><i>Vibrio parahaemolyticus</i></p> <p><i>Vibrio alginolyticus</i></p>
AEROMONADACEAE	<p><i>Aeromonas hydrophila</i>*</p> <p><i>Aeromonas sobria</i></p>	

* hydrolyses ONPG due to an enzyme other than β -galactosidase in the strictest sense of the term, i.e. can be induced by IPTG.

PRECAUTIONS

Respect of Good Laboratory Practice.

BIBLIOGRAPHY

- **RICHARD Cl. (1978):** Annales Biologiques Clin., **36:** 407-424.
- **LE MINOR L., COYNAULT C., GUISON N. (1977):** Annales de l'Institut Pasteur, **128 B:** 35-43.
- **POINDRON P., BONLOUIS C. (1974):** Méd. Mal. Infect., **4:** 23-28.
- **LE MINOR L., BEN HAMIDA F. (1962):** Annales de l'Institut Pasteur, **102:** 267-277.
- **MOLLARET H., LE MINOR L. (1962):** Annales de l'Institut Pasteur, **102:** 649-652.