

## PCB/Agar (Potato-Carrot-Bile)

355-5834

### DEFINITION

Medium used for the identification of yeasts of the genus *Candida*, and in particular *Candida albicans*.

### PRINCIPLE

Nutrient substances favor the growth of yeasts of the genus *Candida*. Due the presence of ox-bile, this medium inhibits bacteria and certain yeasts.

### PRESENTATION

#### Ready-to-use

8 ml x 25 tubes

code 355-5834

### STORAGE

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

### THEORETICAL FORMULA

Carrots, peeled and grated	20 g
Potatoes, peeled and grated	20 g
Bovine bile	150 g
Agar	25 g
Distilled water	1,000 ml
Final pH (25°C) = 6.8 ± 0.2	

### OTHER PRODUCTS REQUIRED (NOT SUPPLIED)

- Distilled water
- **Sabouraud Glucose Agar** (code 356-4494)  
500 g

### EQUIPMENT REQUIRED (NOT SUPPLIED) (non-exhaustive)

- Scales
- Sterile weighing bags
- Grinder
- Sterile pipettes (code 355-0751) or inoculating loops
- Thermostatically-controlled incubator or incubation room, precise to ± 1°C
- All usual laboratory equipment

### PROTOCOL

#### • Inoculation and incubation

Using a Sabouraud glucose agar (code 356-4494) incubated for 24 to 48 hours at 37°C, collect a presumed *Candida* colony and sub-culture it on PCB medium by making several streaks in the bottom of the tube, then a terminal streak at a shallow depth. Incubate at 27°C ± 1°C for 24-48 hours.

### READING AND INTERPRETATION

Use a microscope to detect chlamydo-spores of *Candida albicans* in the filamented zones in which have developed.

If none are present, sub-culture once again on PCB medium and incubate at 27°C for 48 hours.

If still negative, determination of *Candida* species can be identified by using the AUXACOLOR scatter technique and sugar fermentation.

### PRECAUTIONS

- The time lapse between the end of preparation of the stock solution (or the 10<sup>-1</sup> dilution in the case of a solid product) and the moment when the dilutions come into contact with the culture medium must not exceed 15 minutes.
- Comply with Good Laboratory Practice.

### PERFORMANCES / QUALITY CONTROL OF THE TEST

The growth performances of the media are verified with the following strains:

STRAINS	Result after 48 – 72 h at room temperature (*)
<i>Candida albicans</i> ATCC 2091	Positive culture and production of chlamydo-spores
<i>Candida krusei</i>	Positive culture, no chlamydo-spore

(\*) optimum temperature: 27 °C

### 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

PCB (Agar) / *Candida albicans* / Identification / Medium.

**BIBLIOGRAPHY**

- **DROUHET E. - VIEU-COUTEAU M. (December 1957):** Biologie des infections à *Candida*. Diagnostic de laboratoire, Sem. Hôp. Paris.
- **PAVLATOU M. - MARCELOU U. (1956):** Milieu favorisant la formation de chlamydospores de *Candida albicans*. Annales de l'Institut Pasteur 91: 410.