

Blastesis/Agar

355-5844

DEFINITION

Blastesis medium allows the production of "germ tubes" characteristic of *Candida albicans* (98% of strains).

N.B.: *C. stellatoïdea* can also possess this characteristic.

PRESENTATION

Ready-to-use

1.5 ml x 25 tubes

code 355-5844

STORAGE

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

THEORETICAL FORMULA

Synthetic fluid medium.
Final pH (25°C) = 6.8 ± 0.1

PROTOCOL

Surface inoculate the Blastesis medium.
The inoculum must be sufficient to generate very slight turbidity in the medium.
The germ tubes are observed through a microscope, between slide and coverslip, after 3 to 4 hours incubation at 37°C (± 1°C).

PRECAUTIONS

- Take care not to confuse germination with onset of filamentation.
- Comply with Good Laboratory Practice.

PERFORMANCES / QUALITY CONTROL OF THE TEST

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

STRAINS	Production of germ tubes after 3 - 4 h incubation at 37°C	
	T0	T + 3h
<i>Candida albicans</i> ATCC 2091	-	+
<i>Candida albicans</i> ATCC 10231	-	+
<i>Candida tropicalis</i> ATCC 750	-	-

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

Blastesis / *Candida* / Water / Food products / Medium.