

It's Time to *B. cereus*

Introducing Bio-Rad's newest culture media, **RAPID'*B.cereus***

- **Unique** — allows enumeration of all pathogenic *B. cereus* species, including *Bacillus cytotoxicus*
- **Quick turnaround** — produces results in under 24 ± 3 hours with no confirmation required
- **Easy interpretation** — facilitates clear visualization of red colonies surrounded by an opaque halo on a blue background
- **Flexible** — enables surface inoculation and pour plate protocols with storage up to 72 hours at 2–8°C prior to reading

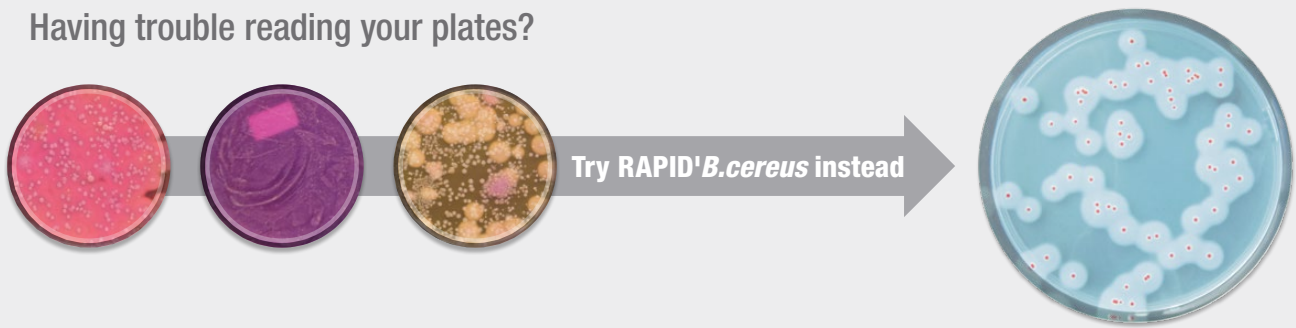
Visit [bio-rad.com/rapidmedia](https://www.bio-rad.com/rapidmedia) for more information and to see how our products can be incorporated into food testing workflows.



BRD 07/26–03/19
ALTERNATIVE ANALYTICAL
METHODS FOR AGRIBUSINESS
Certified by AFNOR Certification
nf-validation.afnor.org/en



Having trouble reading your plates?



Detect all pathogenic *B. cereus* species

RAPID'*B.cereus* is the first chromogenic medium allowing the enumeration of all pathogenic *B. cereus* species, including *Bacillus cytotoxicus*

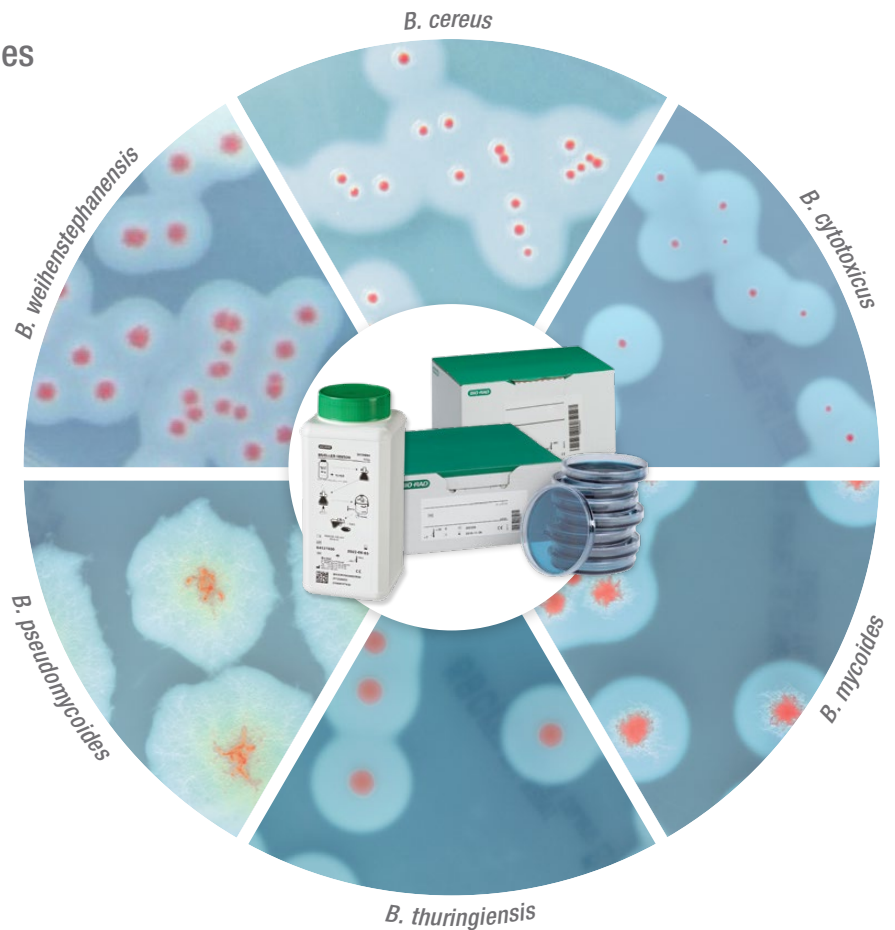
Complete and versatile solution

- Ready-to-use plates
- Dehydrated + supplements
- Surface inoculation
- Pour plate method

Ordering Information

Catalog #	Description
12007304	Ready to use , 90 mm x 20 plates
12007305	Dehydrated base , 500 g
12007306	Supplement 1 , 1 box, 10 vials
12007307	Supplement 2 , 1 box, 10 vials

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

