



RAPID'*P.aeruginosa* Agar for Water Testing

Choose the easy way for your auto-control monitoring
of *Pseudomonas aeruginosa* in bottled water in less than 24 hours



▪ **Easier than conventional methods**

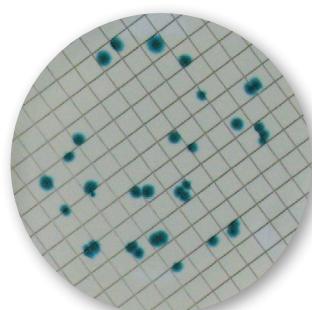
Direct enumeration of *Pseudomonas aeruginosa* in 22-30 hours
No confirmation required

▪ **High enzymatic specificity** delivering highly contrasted
and very specific colony coloration for quick and easy reading

▪ **Validations**

Certified NF VALIDATION as an alternative method
to the NF EN ISO 16266 standard

For more informations: www.foodscience.bio-rad.com



RAPID'*P.aeruginosa* Agar for Water Testing

Principle

RAPID'P.aeruginosa is a selective chromogenic culture medium for the direct enumeration (without confirmation) of *Pseudomonas aeruginosa* using the membrane filtration method.

It is particularly suitable for the testing of water for human consumption, in particular bottled water and untreated water such as water from wells, lakes and streams wherein the commensal water flora is abundant and for which the standardized medium is non-specific for *P. aeruginosa*.

The principle of RAPID'P.aeruginosa medium is based on the detection of an enzymatic activity typical of *P. aeruginosa*.

Under its action, a specific chromogenic substrate is cleaved leading to the formation of a blue to blue-green precipitate on the *Pseudomonas aeruginosa* colonies.

The selective mixture makes it possible to inhibit the majority of interfering flora, very important in water, in particular *Pseudomonas* non *aeruginosa*. Other microorganisms may however show growth.

Their colonies appear transparent or pigmented yellow-green and are easily distinguishable from those of *Pseudomonas aeruginosa*.

Key Benefits

- High specificity for a direct enumeration of *Pseudomonas aeruginosa*
- Complete results within 22-30 hrs
- Reliable results compared to the CN Agar method
- Good colony coloration contrast for rapid assessment
- No confirmation step required

Ordering Information

Catalog # Description

RAPID'P.aeruginosa Agar for Water Testing

Ready to use

356-3984 55 mm x 20 dishes

Dehydrated Base

356-4900 500 g

Detection and enumeration of *Pseudomonas aeruginosa* by membrane-filtration

Certified NF VALIDATION as an alter native method according to the NF EN ISO 16266 standard



Water sample filtration



Membrane

⌚ 22-30 hrs
36°C



RAPID'P.aeruginosa

Enumeration of typical colonies
Blue to blue-green colonies



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 32 3 710 53 00 Brazil 55 31 3689 6600 Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Malaysia 60 3 2117 5260 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 64 9 415 2280 Norway 47 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3170 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 Thailand 66 2 6518311 United Kingdom 020 8328 2000