



## RAPID' *Sakazakii* Agar

**RAPID' *Sakazakii* is a medium for the detection of *Cronobacter* spp.**

- **Easy-to-read chromogenic reaction**

Blue to blue-green colonies

- **Cost-effective short protocol**

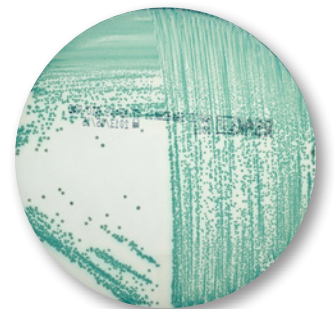
No selective enrichment in mLST + Vancomycine

Negative results in less than 42 hours

- **Validation**

Certified NF VALIDATION according to the ISO 16140 standard for the detection of *Cronobacter* spp. in infant milk powders

For more informations: [www.foodscience.bio-rad.com](http://www.foodscience.bio-rad.com)



RAPID' *Sakazakii* Agar

**The Difference is Clear.** RAPID'Chromogenic Range for Food and Environmental Testing

**BIO-RAD**

## Principle

RAPID'Sakazakii is a selective chromogenic medium used for the detection of *Cronobacter* spp. (formerly *Enterobacter sakazakii*). The principle of the medium is based on the demonstration of an enzymatic activity characteristic of *Cronobacter* spp. Under its action, the chromogenic substrate 5-bromo-4-chloro-3-indolyl- $\alpha$ -D-glucopyranoside is hydrolyzed causing blue to blue-green color for *Cronobacter* spp. colonies. The incubation temperature set at 44°C, sodium deoxycholate and crystal violet inhibit the growth of most of the associated microflora.

To test infant milk powders (very low interfering flora contamination) a validated short protocol can be used to detect *Cronobacter sakazakii* in 24 hours after 24 hours of enrichment in Buffered Peptone Water.

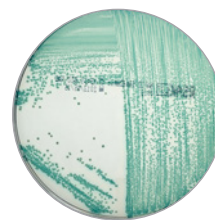
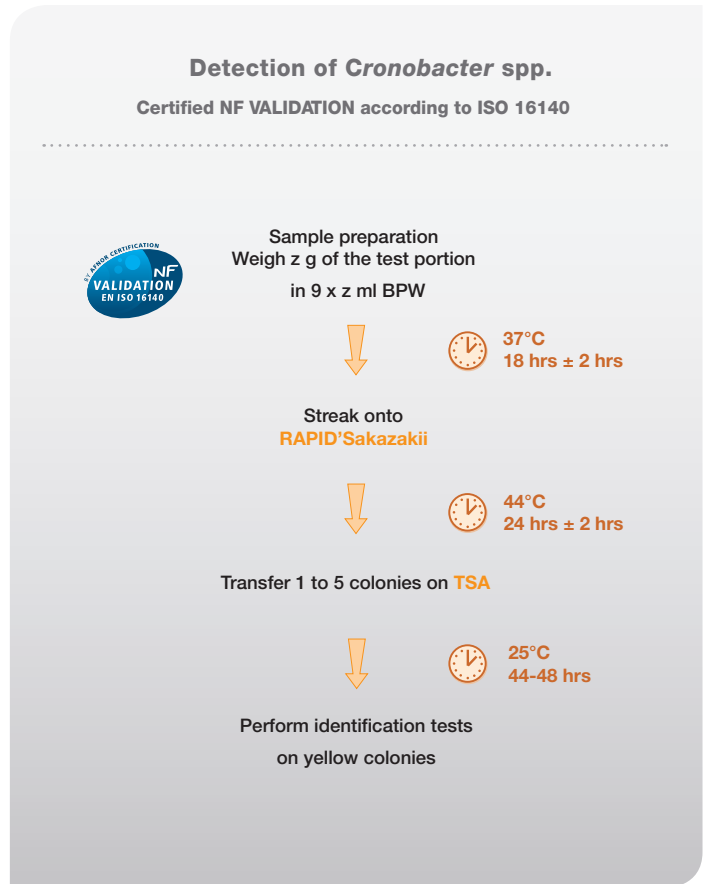
RAPID'Sakazakii chromogenic medium is formulated to be used in ISO 22964, FDA and USDA standard.

## Key Benefits

- Economical: 1 broth, 1 plate
- Fast protocol: negative results in less than 42 hours
- Reliable results: certified NF VALIDATION
- Flexibility: suitable for ISO/TS 22964:2006 standard

## Ordering Information

Catalog #	Description
<b>Buffered Peptone Water/Broth</b>	
<b>Ready to use</b>	
355-4179	225 ml x 6 bottles
355-5795	3 l x 4 bags
355-5790	5 l x 2 bags
<b>Dehydrated</b>	
356-4684	500 g
<b>RAPID'Sakazakii/Agar</b>	
<b>Ready to use</b>	
356-3971	90 mm x 20 dishes
<b>Dehydrated</b>	
356-4976	500 g
<b>Trypto-Casein-Soy (TSA)/Agar</b>	
<b>Ready to use</b>	
356-3884	90 mm x 20 dishes
355-5947	200 ml x 6 bottles
355-5944	8 ml x 25 inclined tubes
<b>Dehydrated</b>	
356-4554	500 g
<b>iQ-Check™Cronobacter/Kit</b>	
357-8137	96 tests



RAPID'Sakazakii



*Cronobacter* spp.



Other  
*Enterobacteriaceae*



**Bio-Rad**  
**Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 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 23 38 41 30 Poland 48 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