

Oxoid™ Cronobacter PreciS™ Method

Publication Number MAN0025440 Revision C00



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Summary

Cronobacter (formerly *Enterobacter sakazakii*) are Gram-negative rod-shaped members of the Family *Enterobacteriaceae* that have been implicated in outbreaks of disease in premature infants (neonates), causing sepsis, meningitis, and necrotizing enterocolitis. Neurological damage can be permanent, and the death rate is reported to be as high as 40–80%.

Cronobacter spp. have been isolated at low levels from powdered infant formulae, and the organisms' high tolerance to desiccation provides a competitive advantage in the dry environments of milk powder factories, increasing the risk of product contamination.

The protocol of the primary enrichment is not clearly described in the ISO 22964:2017 standard for the detection of *Cronobacter* spp. For instance, is there a need to supplement the primary enrichment with antibiotics for infant formula with probiotics or with α -amylase for infant cereals with starch as described in the ISO 6887-4:2017? Additionally, a selective sub-culture is required, extending the time-to-result for 24 hours.

The Oxoid™ Cronobacter PreciS™ Method offers an alternative to the ISO 22964:2017 workflow by requiring only one single enrichment prior to streaking on Oxoid™ Brilliance™ CCI Agar Base (Brilliance™ CCI). This simplified and fast workflow has been validated according to the ISO 16140-2:2016 standard by MicroVal™.

Intended use

Oxoid™ Brilliance™ CCI Agar Base (Brilliance™ CCI) is an optimization of the original Brilliance™ Enterobacter Sakazakii Agar (DFI), developed by Iversen et al. in 2004. The concentration of the X- α -glucoside chromogen was increased to improve the blue-green coloration of *Cronobacter* strains with weak glucosidase activity. Sodium thiosulphate and ammonium iron (III) citrate are included to differentiate hydrogen sulfide-producing *Enterobacteriaceae*, which are colored grey/brown, and sodium deoxycholate is added to inhibit the growth of Gram positive competitors.

Oxoid™ Cronobacter PreciS™ Method offers a fast screening of the samples after an overnight enrichment and a selective plating onto Oxoid™ Brilliance™ CCI Agar Base (Brilliance™ CCI).

This workflow is meant for use in laboratories undertaking microbiological analysis.

Method

The Oxoid™ Cronobacter PreciS™ Method combines the benefits of the Oxoid™ Buffered Peptone Water (ISO) (BPW), Oxoid™ Brilliance™ CCI Agar Base (Brilliance™ CCI), and Oxoid™ Microbact GNB 24E Kit.

This method reduces the time to result over conventional culture methods. The prescribed Oxoid™ BPW is highly nutritious, enabling the fast recovery and growth of *Cronobacter*, including stressed cells.

Oxoid™ Brilliance™ CCI improves recovery of *Cronobacter* by reducing background flora. Chromogens aid easy identification and differentiation by producing brightly colored colonies.

The Oxoid™ Microbact GNB 24E Kit or the SureTect™ Cronobacter species PCR Assay ensure a rapid and simple confirmation for isolated characteristic colonies from the Oxoid™ Brilliance™ CCI plate.

Depending on the legislation territory, it is also possible to use any appropriate EN ISO 16140-6:2019 or Official Method of analysis of AOAC International (AOAC-OMA) validated confirmation method or any appropriate reference confirmation procedure.

Media composition

The following compositions are for typical formulae. Adjustments might be required to meet performance standards.

Table 1 Oxoid™ Buffered Peptone Water (ISO) (Cat. No. [CM1049](#), [CM1211](#), or equivalent)

Reagents	Concentration
Peptone	10 g/L
Sodium Chloride	5 g/L
Disodium Hydrogen Phosphate (anhydrous)	3.5 g/L
Potassium Dihydrogen Phosphate	1.5 g/L
pH 7.0±0.2 at 25°C	

Table 2 Oxoid™ Brilliance™ CCI Base (Cat. No. [CM1122B](#))

Reagents	Concentration
Yeast Extract	3.0 g/L
Sodium Chloride	5.0 g/L
Tryptone	7.0 g/L
Sodium Thiosulphate	1.0 g/L
Sodium Deoxycholate	0.25 g/L
Ferric Ammonium Citrate	1.0 g/L
5-Bromo-4-chloro-3-indolyl α-D-glucopyranoside	0.15 g/L

Table 2 Oxoid Brilliance CCI Base (Cat. No. CM1122B) (continued)

Reagents	Concentration
Agar	13.2 g/L
pH 7.3±0.2 at 25°C	

Materials required (not supplied)

- Inoculating loops, swabs, collection containers
- Incubators
- Quality control organisms

Prepare the materials

Note: Ready-to-use Oxoid™ Buffered Peptone Water (ISO) (Cat. No. BO1067S) can be used as well.

Prepare the Oxoid™ Buffered Peptone Water (ISO)

1. Suspend 4.5 g of Oxoid™ Buffered Peptone Water (ISO) in 225 mL of distilled water.
2. Mix well, distribute into final containers, then sterilize by autoclaving at 121°C for 15 minutes.

Prepare the Novobiocin Selective Supplement

Resuspend or suspend the Oxoid™ Novobiocin Selective Supplement or Oxoid™ PreciSBlue Supplement (Cat. No. [SR0181E](#), [SR0249A](#), [SR0259A](#) or equivalent) as directed in Table 3.

Table 3 Prepare the Oxoid™ Novobiocin Selective Supplement or Oxoid™ PreciSBlue Supplement

Supplement			Reconstitution		Volume of supplement added to volume of BPW (ISO) to achieve 6 mg/L			
Cat. No.	Product name	mg/vial	Solvent	mL/vial	225 mL	1000 mL	1350 mL	3375 mL
SR0259A	Oxoid™ PreciSBlue Selective Supplement (Liquid format)	216	N/A	40	0.25 mL	N/A	1.5 mL	1.04 mL
SR0249A	Oxoid™ Novobiocin Selective Supplement (Liquid format)	400	N/A	40	0.14 mL	0.6 mL	N/A	2.0 mL
SR0181E	Oxoid™ Novobiocin Supplement	10	Distilled sterile water	2	0.27 mL	1.2 mL	N/A	4.1 mL

Prepare the Oxoid™ Vancomycin Supplement

Resuspend the Oxoid™ Vancomycin Supplement (Cat. No. [SR0247E](#) or equivalent) as directed. See recommendation in Table 4.

Table 4 Prepare Oxoid™ Vancomycin Supplement

Supplement			Reconstitution		Volume of supplement added to volume of BPW (ISO) to achieve 6 mg/L		
Cat. No.	Product name	mg/vial	Solvent	mL/vial	225 mL	1000 mL	3375 mL
SR0247E	Oxoid™ Vancomycin Supplement	5	Distilled sterile water	2	0.54 mL	2.4 mL	8.1 mL

Prepare the Oxoid™ *Brilliance™* CCI Agar Base (*Brilliance™* CCI)

1. Suspend 30.6 g of Oxoid™ *Brilliance™* CCI Agar Base (*Brilliance™* CCI) in 1 L of distilled water.
2. Bring to a boil to dissolve completely.
3. Sterilize by autoclaving at 121°C for 15 minutes.
4. Cool to approximately 50°C.
5. Mix well and pour into sterile Petri dishes.

Isolate Cronobacter from infant formula, infant cereals, and related ingredients and from environmental samples

Method certified EN ISO 16140-2:2016 by MicroVal™

Comply with Good Laboratory Practices (refer to EN ISO 7218:2007 standard).

1. Homogenize the sample for 30–60 seconds with a homogenizer.
2. Enrich the samples as follows:

Matrices	Media	Incubation
10 g infant formula with or without probiotics ^[1]	1:10 ratio of sample to media. For example, add up to 10 g of sample to 90 mL of Oxoid™ Buffered Peptone Water (ISO)	34°C to 38°C for 16–22 hours
Up to 375 g infant formula with and without probiotics, infant cereals with and without probiotics, and related ingredients ^[1]	1:10 ratio of sample to media. For example, add up to 375 g of sample to 3,375 mL of Oxoid™ Buffered Peptone Water (ISO) supplemented with 6 mg/L of vancomycin solution	34°C to 38°C for 20–26 hours
	1:6 ratio of sample to media. For example, 375 g of sample and 1,875 mL of pre-warmed Buffered Peptone Water (ISO) (supplemented with 6 mg/L novobiocin/PrecisBlue for probiotics only)	34°C to 38°C for 18–24 hours
Environmental samples	1:10 ratio of sample to media <ul style="list-style-type: none"> • Add up to 25 g of sample to 225 mL of Oxoid™ Buffered Peptone Water (ISO) supplemented with 6 mg/L of vancomycin solution • Add one swab to 10 mL of Oxoid™ Buffered Peptone Water (ISO) supplemented with 6 mg/L of vancomycin solution • Add one sponge to 100 mL of Oxoid™ Buffered Peptone Water (ISO) supplemented with 6 mg/L of vancomycin solution 	34°C to 38°C for 18–24 hours
	1:10 ratio of sample to media <ul style="list-style-type: none"> • Add up to 25 g of sample to 225 mL of Oxoid™ Buffered Peptone Water (ISO) • Add one swab to 10 mL of Oxoid™ Buffered Peptone Water (ISO) • Add one sponge to 100 mL of Oxoid™ Buffered Peptone Water (ISO) 	34°C to 38°C for 20–26 hours

^[1] The enrichment conditions are harmonized with Oxoid™ Salmonella PreciS™, SureTect™ Salmonella species PCR Assay, and SureTect™ Cronobacter species PCR Assay, with no addition of α-amylase required.

3. Gently agitate the bag. Then, using a microbiological loop, inoculate a 10 µL loopful of the broth onto a plate of Oxoid™ *Brilliance*™ CCI Agar Base (*Brilliance*™ CCI) using a diminishing sweep technique to produce single colonies.

Note: It is possible to store the enrichment broth after incubation for 72 hours at 5°C±3°C before streaking onto *Brilliance*™ CCI except when testing 10 g infant formula with or without probiotics enriched in non-supplemented Oxoid™ Buffered Peptone Water.

4. Incubate the plates at 37±1°C or 41.5±1°C for 24±2 hours.

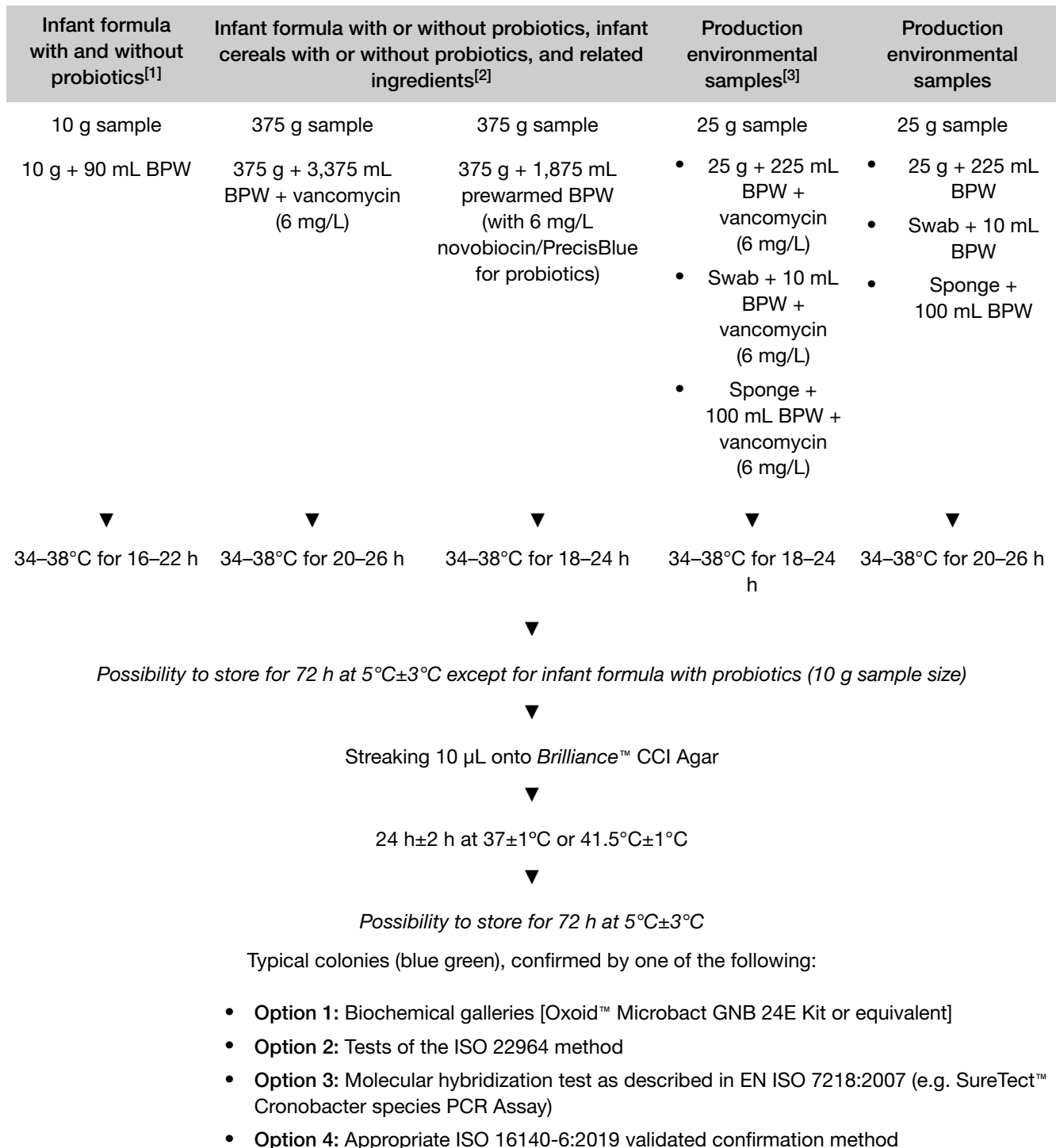
Note: It is possible to store the plates for 72 hours at 5°C±3°C before confirmation.

Cronobacter colonies grow as 1–3 mm blue/green colonies, and non-target organisms are either inhibited or grow as straw or white colonies.

Blue colonies are presumptive positive for *Cronobacter* (see Figure 1).

See Confirm positive results on page 9 to confirm the observed characteristic colonies.

Oxoid™ Cronobacter Precis™ Method Workflow



^[1] No addition of α-amylase required.

^[2] No addition of α-amylase required. It is possible to test smaller sample sizes by adjusting the volume of BPW + vancomycin (6 mg/L).

^[3] For sampling after cleaning process, premoisten:

- 1 swab + 1 mL broth universal neutralizing (+ 9 mL BPW)
- 1 sponge + 10 mL broth universal neutralizing (+ 90 mL BPW)

- 1 wipe + BPW + 10% neutralizing agent (+ 225 mL BPW)

Example results

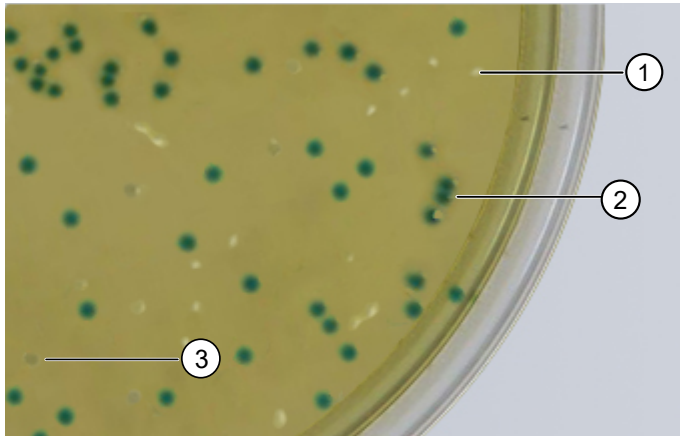


Figure 1 Example results—Mixed culture

- ① *E. coli* colony
- ② *Cronobacter spp.* colony
- ③ *Enterobacter cloacae* colony

Confirm positive results



Confirmation is performed from isolated characteristic colonies on Oxoid™ Brilliance™ CCI Agar Base (Brilliance™ CCI) and running one of these options:

- Option 1: Biochemical galleries [Oxoid™ Microbact GNB 24E Kit (Cat. No. [MB1131A](#)) or equivalent].
- Option 2: EN ISO 22964:2017 confirmation procedure.
- Option 3: Any appropriate EN ISO 16140-6:2019 validated method.
- Option 4: Molecular hybridization as described in EN ISO 7218:2007 using for instance the SureTect™ Cronobacter species PCR Assay (Cat. No. [PT1060A](#) or [A56845](#)) that is validated according to the ISO 16140-2:2016 standard (NF Validation certificate UNI 3/11-12/15).

In the event of discordant results (positive with the Oxoid™ Cronobacter Precis™ Method or non-confirmed by one of the means described above), the laboratory must follow the necessary steps to ensure the validity of the result obtained.

Performance validation

Table 5 MicroVal™ certification of the method

Certification	Scope and expiration
 <p>MICROVAL®  NEN</p> <p>MicroVal™ certificate 2020LR93</p> <p>http://www.microval.org</p>	<p>The MicroVal™ certificate can be obtained from our technical support team.</p> <ul style="list-style-type: none">• email: microbiology.techsupport.uk@thermofisher.com• telephone: +44 (0)1256 694238• or from the MicroVal™ website: http://www.microval.org <p>For more information about the validity of the MicroVal™ certification, see the certificate 2020LR93 available at http://www.microval.org or obtain from our technical support team.</p>

References

- Iversen C., Druggan P., Schumacher S., Lehner A., Feer C., Gschwend K., Joosten H., Stephan R., 2008. Development of a novel screening method for the isolation of *Cronobacter* spp. (*Enterobacter sakazakii*). *Appl. Environ. Microbiol.* 74:2550–2552.
- Mullane N. R., Iversen C., Healy B., Walsh C., Whyte P., Wall P. G., et al, 2007. *Enterobacter sakazakii* an emerging bacterial pathogen with implications for infant health. *Minerva Pediatr.* 59:137–148.
- Lai K.K., 2001 *Enterobacter sakazakii* infections among neonates, infants, children, and adults: case reports and a review of the literature. *Medicine (Baltimore)* 80:113–122.
- EN ISO 22964:2017. Microbiology of the food chain – Horizontal method for the detection of *Cronobacter* spp.
- EN ISO 7218:2007. Microbiology of food and animal feeding stuffs – General requirements and guidance for microbiological examinations.
- EN ISO 16140-2:2016. Microbiology of food and animal feed – Method validation – Part 2: Protocol for the validation of alternative (proprietary) methods against a reference method.
- EN ISO 16140-6:2016. Microbiology of food and animal feed – Method validation – Part 6: Protocol for the validation of alternative (proprietary) methods for microbiological confirmation and typing procedures.

Revision history: MAN0025440 C00 (English)

Revision	Date	Description
C00	8 April 2024	A new Novobiocin selective supplement was added, Oxoid™ PrecisBlue Selective Supplement.
B.0	14 December 2023	<ul style="list-style-type: none">• Matrices were added for infant formula with and without probiotics, infant cereals with and without probiotics, and related ingredients.• Matrices were added for environmental samples in Oxoid™ Buffered Peptone Water (ISO) not supplemented with vancomycin solution.• Instructions were added for how to prepare Novobiocin.
A.0	28 September 2021	New document for MicroVal™ validation study of Oxoid™ Cronobacter Precis™ Method.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Limited Use Label License: This product or portions hereof are subject to proprietary rights of Epoch Biosciences, Inc. and are made and sold under license from Epoch under the patents and patent applications as may be designated by Epoch from time to time set forth, including one or more of the following: U.S. Patent Nos. 6,699,975, 7,718,374, 7,348,146 and 7,671,218, and related U.S. and foreign patents and pending applications. A license must be obtained directly from Epoch with respect to any other proposed use of this product.

Buyer agrees and understands it is not licensed to (a) provide commercial data or databases to or on behalf of any third party that relate to use of the product; or (b) use the product or data there from in a clinical diagnostic setting where data from an individual's sample is given to such individual or their caregiver. Neither Epoch Biosciences nor its licensors grants any other licenses, express or implied, to permit the manufacture, use, sale, or importation of the product, any other nucleic acids, or any data generated from use of the product.

Trademarks: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. MicroVal is a trademark of MicroVal.

©2021-2024 Thermo Fisher Scientific Inc. All rights reserved.