



PIF Supplement Workflow Benefits and Performance Data

BIO-RAD

Bio-Rad Laboratories

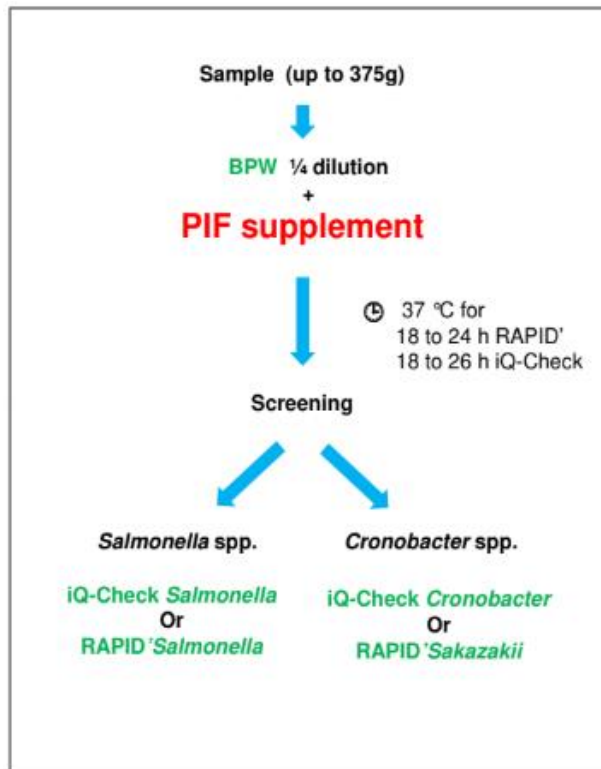
Food for Thought Newsletter, June 2020



Workflow Benefits

BIO-RAD

A Fully Harmonized Solution



Harmonized

Infant formula without probiotics
Infant formula with probiotics
Ingredients



Harmonized

Salmonella
Cronobacter



Harmonized

iQ-Check Kits
RAPID' media

Benefits of 1:4 Dilution

Make your technicians happier with 1:4 dilution

375 g

1:10



3.75 kg / bag



375 g

1:6



2.25 kg / bag



375 g

1:4



1.5 kg / bag



Workflow Benefits

A small 1:4 dilution factor for big benefits



Summary

PERFORMANCE

Very low level of detection, even in
challenging conditions:
Large sample size
Stressed bacteria

EASE OF USE

Large sample size suitable method
No additional materials / technicians
Co-enrichment *Salmonella* & *Cronobacter*

REASONABLE COST

Reagent cost
Consumable & material cost
Labor cost



Performance Data

AFNOR Validation Data
RAPID' *Sakazakii*

BIO-RAD

RAPID' *Sakazakii* AFNOR Performance Data

	Category 1: Infant formula and infant cereals with and without probiotics including ingredients (30 g sample size)	Category 2: Infant formula and infant cereals with and without probiotics including ingredients (375 g sample size)
Enrichment	30 g + 270 mL BPW* 18 h \pm 2 h at 37°C \pm 1°C	375 g + 1125 mL Pre-warmed (37°C) BPW with PIF supplement* 18 to 24 h at 37°C \pm 1°C
Streaking	10 μ L onto RAPID' <i>Sakazakii</i> Incubation 24 h \pm 2 h at 44°C \pm 1°C	
Confirmation	1 to 5 colonies by the tests described in the ISO method (biochemical tests) with or without purification step on TSA or by PCR (iQ Check <i>Cronobacter</i>) on isolated colony	
Enrichment broth storage	48 h at 5°C \pm 3°C	72 h at 5°C \pm 3°C
Plate storage	72 h at 5°C \pm 3°C	

RAPID[®] Sakazakii AFNOR Performance Data

Sensitivity Study

Category	Type	SE alt %	SE ref %	RT %	FP %
Infant formula and infant cereals with and without probiotics including ingredients (30 g sample size)	Infant formula and infant cereals without probiotics	100,0	100,0	100,0	0,0
	Infant formula and infant cereals with probiotics	100,0	100,0	100,0	0,0
	Ingredients (Maltodextrin, NFDM, whey...)	85,7	100,0	90,0	0,0
	Total	94,6	100,0	96,7	0,0
Infant formula and infant cereals with and without probiotics and ingredients (375 g sample size)	Infant formula and infant cereals without probiotics	100,0	84,6	90,9	0,0
	Infant formula and infant cereals with probiotics	100,0	81,8	90,9	0,0
	Ingredients (Maltodextrin, NFDM, whey...)	100,0	50,0	77,3	0,0
	Total	100,0	73,5	86,4	0,0
Total		97,2	87,3	91,3	0,0

RLOD Study

Category	Name	Study design	AL	RLOD
1	Infant cereals with probiotics / <u><i>Cronobacter sakazakii</i></u> QL 11007.9	Unpaired	2,5	0,843
2	Infant formula with probiotics / <u><i>Cronobacter sakazakii</i></u> Ad1418			0,868
Combined				0,954

Better sensitivity than the reference method

No false positive results

RAPID'Sakazakii AFNOR Performance Data

AFNOR certificate (available)



Contacts :

Méline TERRO
Ingénierie certification
meline.terro@afnor.org
Tél : +33 (0)1 41 62 62 39

Réf.: MET/MET/NF102/Clients/BIO-RAD/
Avis BT_RAPID'Sakazakii_2020-04-02 (R2).docx

BIO-RAD
M. Yannick BICHOT
3 boulevard Raymond Poincaré
92430 MARNES-LA-COQUETTE

Objet : Marque NF VALIDATION

La Plaine Saint-Denis, le 2 avril 2020

Monsieur,

Comme suite à l'avis positif exprimé le 2 avril 2020 par le Bureau Technique de la marque NF VALIDATION (NF102), dans son application à l'agroalimentaire, j'ai l'honneur de vous annoncer que le **droit d'usage** de la marque NF VALIDATION est reconduit pour la méthode alternative suivante :

RAPID'Sakazakii

Certifiée sous le N° BRD 07/22-05/12, avec pour fin de validité le 10-Mai-2024

La méthode alternative est validée par comparaison à la méthode de référence NF EN ISO 22964 (2017) et selon le protocole de validation NF EN ISO 16140-2 (2016) pour la détection des *Cronobacter* spp. dans le domaine d'application suivant : « poudres de lait infantile et céréales avec ou sans probiotiques infantiles incluant leurs ingrédients (avec prise d'essai de 30g) », « poudres de lait infantile et céréales infantiles avec ou sans probiotiques incluant leurs ingrédients (avec prise d'essai de 375g) ».

Un courrier complet de conclusions mentionnant d'éventuelles réserves prononcées par le Bureau Technique, vous sera prochainement adressé. Le cas échéant, celles-ci devront être prises en compte sans délai.

Je vous prie d'agréer, Monsieur, mes salutations distinguées.

Directeur Général
Franck LEBEUGLE

11 rue Francis de Pressensé - 93571 La Plaine Saint-Denis Cedex - France - T : +33 (0)1 41 62 80 00 - F : +33 (0)1 49 17 90 00
SAS au capital de 18 167 000 € - 479 076 002 RCS Bobigny - www.afnor.org



CERTIFY THE ANALYTICAL PERFORMANCES OF TEST KITS

Validation Summary Report (soon available)



NF VALIDATION
Validation of alternative analytical methods
Application in food microbiology

Summary Report

Validation study according to the EN ISO 16140-2:2016

RAPID'Sakazakii

(Certificate number: BRD 07/22 - 05/12)

for the detection of *Cronobacter* spp. in infant formula and infant cereals with or without probiotics including ingredients (30 g and 375 g sample sizes)

Qualitative method

Expert Laboratory:	ADRIA Développement ZA Creach Gwen 29196 Quimper Cedex (France)
For:	BIO-RAD 3 boulevard Raymond Poincaré 92430 Marnes la Coquette (France)

This report consists of 75 pages, including 7 appendices.
Only copies including the totality of this report are authorised.
Competencies of the laboratory are certified by COFRAC accreditation for the analyses marked with the symbol*.

Version 0
30 April 2020

ADRIA DEVELOPPEMENT • Creach Gwen • F. 29196 QUIMPER Cedex • Tél. (33) 02 98 10 18 18 • Fax (33) 02 98 10 18 99
adria.developpement@adria.tm.fr • <http://www.adria.tm.fr>
ASSOCIATION LOI DE 1905 - N° SIRET 305 964 271 00086 - N° EXISTENCE 53200006329 - N° TVA FR48309964271



Performance Data

AFNOR Validation Data
iQ-Check Cronobacter

BIO-RAD

iQ-Check *Cronobacter* AFNOR Performance Data

Step	Category 1: Infant formula and infant cereals with and without probiotics including ingredients	Category 2: Infant formula and infant cereals with and without probiotics including ingredients	Category 3: Environmental samples
Enrichment	30 g + 270 mL BPW + 10 µg/mL vancomycin 20h ± 2h at 37°C ± 1°C Subculture: 1mL enriched BPW + 9mL BPW 4h ± 1h at 37°C ± 1°C	375 g + 1125 mL Pre- warmed (37°C) BPW with PIF supplement* 18 h to 26 h at 37°C ± 1°C	30g + 270 mL BPW or Sampling device + 10 mL (Swab), + 100 mL (sponge) + 225 mL (wipe) 18h +/- 2h at 37°C
FDRS protocol**	Applied on	100 µl of decanted enriched broth.	
DNA extraction	Easy extraction protocol or 100µL	using Deepwell plates on FDRS treated enriched sample	100µL enriched sample
PCR	On 5 µL DNA extract with the CFX Manager software IDE	using the CFX96 Real-time Thermocycler version 3.0	
Confirmation of PCR tests	Streaking 10 µL of last enriched sample (BPW) onto RAPID [®] Sakazakii Incubation 24 h ± 2 h at 44°C ± 1°C	Streaking 10 µL onto RAPID [®] Sakazakii Incubation 24h ± 2h at 44°C ± 1°C	Subculture of BPW in mLST (0.1 mL + 10 mL) for 24 h ± 2 h at 44°C ± 1°C Streaking 10 µL onto RAPID [®] Sakazakii Incubation 24h ± 2h at 44°C ± 1°C
Confirmation of typical colonies	1 to 5 colonies by the tests described in the ISO method (Oxidase and biochemical tests) after purification step on non-selective agar plate or without purification step from RAPID [®] Sakazakii (biochemical tests)		
Enrichment broth storage	48 h at 5°C ± 3°C	72 h at 5°C ± 3°C	48 h at 5°C ± 3°C

* Addition of α-amylase for infant cereals (0.1 g/L)

** Optional

- **AFNOR validation extension:**
 - 375 g sample size
 - 1:4 dilution in BPW pre-warmed + PIF Supplement
 - 18–26 hr single enrichment @ 37 ± 1°C
 - Free DNA Removal Solution (FDRS)
 - APF Fast: “Crono Fast”
 - 72 hr storage
- **Matrices**
 - Infant formula and cereals w/wo probiotics and ingredients
 - Environmental samples

iQ-Check *Cronobacter* AFNOR Performance Data

Sensitivity Study

Category	SE Alt %	SE Ref %	RT %	FP %
1 – 30 g regrowing	94.6	100.0	96.7	8.7
2 – 375 g + PIF	100	73.5	86.4	0.0
3 – Environmental	100.0	100.0	100.0	0.0

Sensitivity for the alternative method	$SE_{alt} = \frac{(PA + PD)}{(PA + ND + PD)} \times 100\%$	98.1 %
Sensitivity for the reference method	$SE_{ref} = \frac{(PA + ND)}{(PA + ND + PD)} \times 100\%$	91.4 %
Relative trueness	$RT = \frac{(PA + NA)}{N} \times 100\%$	94.3 %
False positive ratio for the alternative method* FP = PPNA + PPND	$FPR = \frac{(FP)}{NA} \times 100\%$	2.3 %

With $ND = ND + PPND$
 $NA = NA + PPNA$

iQ-Check *Cronobacter* AFNOR Performance Data

RLOD

1 – 30 g regrowing

2 – 375 g + PIF

3 – Environmental

Category	(Strain / matrix) pair	Level of detection at 50% (CFU / sample size) according to Wilrich & Wilrich ¹	
		Reference method	Alternative method
1	Infant cereals with probiotics / <i>Cronobacter sakazakii</i> QL 11007.9	1,5 [0,8;2,9]	1,3 [0,7;2,5]
2	Infant formula with probiotics / <i>Cronobacter sakazakii</i> Ad1418	1,0 [0,6;1,8]	0,9 [0,5;1,6]
3	Process water / <i>Cronobacter sakazakii</i> CCUG 28868	1,5 [0,8;2,6]	1,7 [0,9;3,0]
Combined results		1,3 [0,9;1,8]	1,2 [0,9;1,7]

Category	Name	Study design	AL	RLOD	RLODL	RLODU	b=ln(RLOD)	sd(b)	z-Test statistic	p-value
1	Infant cereals with probiotics / <i>Cronobacter sakazakii</i> QL 11007.9	Unpaired	2,5	0,843	0,297	2,397	-0,17	0,522	0,326	1,256
2	Infant formula with probiotics / <i>Cronobacter sakazakii</i> Ad1418			0,868	0,392	1,926	-0,141	0,398	0,354	1,277
3	Process water / <i>Cronobacter sakazakii</i> CCUG 28868			1,159	0,455	2,953	0,148	0,467	0,316	0,752
Combined				0,954	0,57	1,598	-0,047	0,258	0,183	1,145

iQ-Check *Cronobacter* AFNOR Performance Data

AFNOR certificate (available)



Contacts :

Méline TERRO
Ingénierie certification
melanie.terro@afnor.org
Tél : +33 (0)1 41 62 62 39

Ref: MET/MET/NF102/Clients/BIO-RAD/
Avis BT_iQ-Check Crono_2020-04-02 (R2).docx

BIO-RAD
M. Jean-Philippe TOURNIAIRE
3 boulevard Raymond Poincaré
92430 MARNES-LA-COQUETTE

Objet : Marque NF VALIDATION

La Plaine Saint-Denis, le 2 avril 2020

Monsieur,

Comme suite à l'avis postif exprimé le 2 avril 2020 par le Bureau Technique de la marque NF VALIDATION (NF102), dans son application à l'agroalimentaire, j'ai l'honneur de vous annoncer que le **droit d'usage** de la marque NF VALIDATION est reconduit pour la méthode alternative suivante :

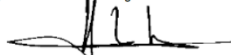
iQ-Check® *Cronobacter* spp.

Certifiée sous le N° BRD 07/23-01/13, avec pour fin de validité le 31-Jan.-2025

La méthode alternative est validée par comparaison à la méthode de référence NF EN ISO 22964 (2017) et selon le protocole de validation NF EN ISO 16140-2 (2016) pour la détection des *Cronobacter* spp. dans les poudres de lait infantile et céréales infantiles avec ou sans probiotiques incluant leurs ingrédients (avec prise d'essai de 30g), poudres de lait infantile et céréales infantiles avec ou sans probiotiques incluant leurs ingrédients (avec prise d'essai de 375g) et les échantillons de l'environnement de production industrielle. Cette reconduction a permis d'étendre le domaine d'application à l'utilisation d'un nouveau protocole d'enrichissement pour les prises d'essais de 375g associé au protocole de lyse Easy, le protocole iQ-Check Free DNA Removal Solution (FDRS) associé au protocole de lyse Easy et le thermoprofile PCR raccourci (APF Fast pour les méthodes iQ-Check) « Crono Fast ».

Un courrier complet de conclusions mentionnant d'éventuelles réserves prononcées par le Bureau Technique, vous sera prochainement adressé. Le cas échéant, celles-ci devront être prises en compte sans délai.

Je vous prie d'agréer, Monsieur, mes salutations distinguées.


Directeur Général
Franck LEBEUGLE

11 rue Francis de Pressensé - 93571 La Plaine Saint-Denis Cedex - France - T: +33 (0)1 41 62 80 00 - F: +33 (0)1 49 17 90 00
SAS au capital de 19 187 000 € - 479 076 002 RCS Bobigny - www.afnor.org



CERTIFY THE ANALYTICAL PERFORMANCES OF TEST KITS

Validation Summary Report (soon available)



NF VALIDATION
Validation of alternative analytical methods
Application in food microbiology

Summary Report

Validation study according to the EN ISO 16140-2:2016

iQ-Check® *Cronobacter* spp. method

(Certificate number: BRD 07/23-01/13)

for the detection of *Cronobacter* spp. in infant formula and infant cereals with or without probiotics including ingredients (30 g and 375 g) and environmental samples

Qualitative method

Expert Laboratory:	ADRIA Développement ZA Creach Gwen 29196 Quimper Cedex (France)
For:	BIO-RAD 3 boulevard Raymond Poincaré 92430 Marnes la Coquette (France)

This report consists of 86 pages, including 7 appendices.
Only copies including the totality of this report are authorized.
Competencies of the laboratory are certified by COFRAC accreditation for the analyses marked with the symbol*.

Version 0
29 April 2020

ADRIA DEVELOPPEMENT • Creach Gwen - F. 29196 QUIMPER Cedex • Tél. (33) 02 98 10 18 18 • Fax (33) 02 98 10 18 39
adria.developpement@adria.fr • <http://www.adria.fr>
ASSOCIATION L29 02 1987 - N° SIRET 339 084 271 9008 - N° COFRAC 520309020 - N° IVA F063396421

Thank You

