



PIF Supplement — A Fully Harmonized Solution

The PIF Supplement has been designed for infant formula, milk powders, cereals, and related ingredients for producers and service laboratories that demand high-performance methods for *Cronobacter*, *Enterobacteriaceae*, and *Salmonella* detection in large sample (up to 375 g) analysis. Our product meets this market segment's strict quality requirements and provides unmatched performance in an easy-to-use format, offering security and peace of mind.


- Single enrichment broth at 1:4 dilution and homogeneous protocol for ease of use
- Product applications for both iQ-Check PCR Detection Kit and RAPID'Chromogenic Media methods
- Validated methods*

* Check with your Bio-Rad representative for validation information.

Visit bio-rad.com/pif for more information and to see how the PIF Supplement can be incorporated into food safety testing workflows.

PIF Supplement Enrichment

BPW + PIF



Up to 375 g:

- Infant formula with or without probiotics
- Milk powders
- Cereals
- Related ingredients

Enrich →

iQ-Check *Cronobacter* spp. Kit







iQ-Check *Enterobacteriaceae* Kit

iQ-Check *Salmonella* II Kit

RAPID' *Sakazakii* Agar

RAPID' *Salmonella* Agar

Save Time, Money, and Resources with PIF Supplement

375 g/1:4 Dilution	375 g/1:6 Dilution	375 g/1:10 Dilution
<p>BPW + PIF</p>  <p>1,125 ml BPW Total enrichment weight = 1.5 kg</p> <p>↓</p> <p>✓ <i>Cronobacter</i> ✓ <i>Enterobacteriaceae</i> ✓ <i>Salmonella</i></p> <p>FULLY harmonized for <i>Cronobacter</i>, <i>Enterobacteriaceae</i>, and <i>Salmonella</i>. Identical conditions and protocols for all three targets.</p> 	<p>Enrichment 1</p>  <p>1,875 ml BPW Total enrichment weight = 2.25 kg</p> <p>↓</p> <p>? <i>Cronobacter</i> ? <i>Enterobacteriaceae</i> ? <i>Salmonella</i></p> <p>NOT FULLY harmonized for <i>Cronobacter</i>, <i>Enterobacteriaceae</i>, and <i>Salmonella</i>. Conditions and protocols need to be verified.</p> 	<p>Enrichment 2</p>  <p>3,375 ml BPW Total enrichment weight = 3.75 kg</p> <p>↓</p> <p>? <i>Cronobacter</i> ? <i>Enterobacteriaceae</i> ? <i>Salmonella</i></p> <p>NOT FULLY harmonized for <i>Cronobacter</i>, <i>Enterobacteriaceae</i>, and <i>Salmonella</i>. Conditions and protocols need to be verified.</p> 
	+66%	+200%
👍	👎	👎

Ordering Information

Catalog #	Description	Catalog #	Description	Catalog #	Description	Catalog #	Description
PIF Supplement		BPW Plus		iQ-Check <i>Cronobacter</i> spp. Kit		RAPID[®] <i>Sakazakii</i> Agar	
12013322	2 g vial	Ready to use		3578137	96 tests	Ready to use	
BPW Standard		3554179	225 ml x 6 bottles	iQ-Check <i>Enterobacteriaceae</i> Kit		3563971	90 mm x 20 dishes
Ready to use		3555795	3 L x 4 bags	12003068	96 tests	Dehydrated	
12013260	5 L x 2 bags	3555790	5 L x 2 bags	iQ-Check <i>Salmonella</i> II Kit		3564976	500 g
Dehydrated		Dehydrated		3578123	96 tests	RAPID[®] <i>Salmonella</i> Agar	
12013259	500 g	3564684	500 g	iQ-Check Free DNA Removal Solution		Ready to use	
12013258	5 kg	3564686	5 kg	3594970	175 mg	3563961	90 mm x 20 dishes
						3563963	90 mm x 120 dishes
						Dehydrated	
						3564705	500 g

Visit bio-rad.com/iqcheck and bio-rad.com/rapidmedia for additional related product information.

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



Bio-Rad
Laboratories, Inc.

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

