

Buffered Peptone Water Plus

Catalog #	Description
3554101	BPW Plus , 100 ml x 6 bottles
3554179	BPW Plus , 225 ml x 6 bottles
3555790	BPW Plus , 5 L x 2 bags
3555795	BPW Plus , 3 L x 4 bags
3564684	BPW Plus , dehydrated, 500 g
3564686	BPW Plus , dehydrated, 5 kg

For laboratory use only.

Intended Use

Buffered Peptone Water (BPW) Plus is used as diluent and nutritive enrichment broth for the resuscitation and growth of a wide variety of microorganisms including *Cronobacter*, *Enterobacteriaceae*, *Escherichia coli*, and *Salmonella* in all food products, animal feed products, and water samples. Notably, it is adapted to demanding strains, stringent growth conditions, and challenging protocols.

Principle

The special selection of peptones allows a very high level of nutritivity to ensure the best recovery. In combination with the phosphate buffer's maintenance of a steady pH and osmotic balance, thanks to the presence of chloride sodium, BPW Plus offers favorable conditions for the growth of a wide range of microorganisms from a variety of matrices.

Theoretical Composition

Peptone 10 g

Sodium chloride 5 g

Potassium dihydrogenphosphate 1.5 g

Disodium hydrogen phosphate 3.5 g

Final pH at 25°C = 7.0 ± 0.2

Shelf Life and Storage

Store dehydrated media at 15–25°C in a carefully sealed package in a dry and dark place.

Store liquid media at 15–25°C.

Required Materials Not Supplied

This is a non-exhaustive list.

Equipment

- All usual laboratory equipment
- Incubators or incubation room
- Scales
- Stirrer/homogenizer
- Vortexer

Precautions

- Respect Good Laboratory Practice (EN ISO 7218). Appropriate protection, such as gloves and lab coats, should be worn when working with potentially infectious live bacteria
- RAPID[®]chromogenic and iQ-Check methods that contain an enrichment in BPW are developed and validated with BPW Plus for optimal performance of the method
- Media that have come in contact with food samples should be considered contaminated and should be disposed of in accordance with local rules and regulations
- For SDS product safety information and certificate of analysis, visit bio-rad.com

Quality Control

Every product manufactured and marketed by Bio-Rad is subject to a quality assurance procedure at all stages, from reception of raw materials through to marketing of the finished products. Each batch of finished product undergoes quality control according to EN ISO 11133 and is marketed only if it satisfies the acceptability criteria. Documentation relative to the production and quality control of each batch is kept on file.

Protocol

Dehydrated BPW Plus Preparation

- Dissolve 20 g of BPW Plus in 1,000 ml of sterile distilled water
- Mix, heating if necessary, until a homogeneous suspension is obtained
- Dispense in appropriate container
- Sterilize in autoclave at $121 \pm 1^\circ\text{C}$ for 15 min

Sample Preparation and Enrichment Protocol

- Dilute sample according to the standard method applicable to the product concerned

References

ISO 11290-2:2017 – Microbiology of the food chain – Horizontal method for the detection and enumeration of *Listeria monocytogenes* and of *Listeria* spp. – Part 2: enumeration method

ISO 19250:2013 – Water quality – Detection of *Salmonella* spp.

ISO 21528-1:2017 – Microbiology of the food chain – Horizontal methods for the detection and enumeration of *Enterobacteriaceae* – Part 1: detection of *Enterobacteriaceae*

ISO 22964:2017 – Microbiology of the food chain – Horizontal method for the detection of *Cronobacter* spp.

ISO 6579-1:2017 – Microbiology of the food chain – Horizontal method for the detection, enumeration and serotyping of *Salmonella* – Part 1: Detection of *Salmonella* spp.

ISO 6887-1:2017 – Microbiology of the food chain – Preparation of test samples, initial suspension and decimal dilutions for microbiological examination – Part 1: General rules for the preparation of the initial suspension and decimal dilutions

ISO 6887-5:2010 – Microbiology of food and animal feeding stuffs – Preparation of test samples, initial suspension and decimal dilutions for microbiological examination – Part 5: specific rules for the preparation of milk and milk products

Revision History

Release date	Document number	Change
06-12-2020	10000128740 Ver A	<ul style="list-style-type: none"> - Major change - Product name change - New document design - Document number change – previous version V6_10/04/2013

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.
All trademarks used herein are the property of their respective owner.