

SOLUS ONE SUPPLEMENT

Insert Number: 44
Issue Number: 1.0
Date of Issue: January 2020
Product Code(s): SALSUPP22.5
SALSUPP112.5



1. PRODUCT INFORMATION

Solus One Supplement is a supplement for the selective enrichment of Gram negative bacteria in Solus One assay testing.

2. SOLUS ONE SUPPLEMENT PACK SIZES

	SALSUPP22.5	SALSUPP112.5
Number of vials	10	10
Volume of 70% Ethanol v:v required per vial	10ml	50ml
“Full strength” (supplement required per L of media)	4.44ml/L	4.44ml/L
“Half strength” (supplement required per L of media)	2.22ml/L	2.22ml/L

3. METHOD FOR RECONSTITUTION

- 3.1. Prepare enrichment media (e.g. BPW) according to manufacturer’s instructions. It is recommended to test the enrichment media to ensure it supports the growth of the target organism according to ISO 11133.
- 3.2. Pre-warm 70% Ethanol v:v to 41.5°C. For reference, a 500ml bottle requires approximately 2 hours in 41.5°C incubator to achieve temperature (from room temperature starting point).
- 3.3. Remove the plastic cap from the supplement vial and discard appropriately.
- 3.4. Using aseptic technique, remove the rubber bung from the vial being careful to avoid any cross contamination.
- 3.5. Aseptically add 70% Ethanol v:v to the vial (see section 2 for volume required in each vial) and allow contents to dissolve into solution.
- 3.6. Once fully dissolved into solution, supplement enrichment media to the level appropriate for the matrix and pathogen to be tested. See individual assay instructions for further information.

4. STORAGE OF RECONSTITUTED SUPPLEMENT

Replace the rubber bung into the vial and store at 2-8°C for up to 2 weeks. Sediment may settle to the bottom of the vial whilst in storage. The efficacy of the supplement is not affected by the presence of this sediment.

5. STORAGE OF SUPPLEMENTED ENRICHMENT MEDIA

A capped container can be stored at 2-8°C for up to 2 weeks. Ensure supplemented enrichment media is brought to the correct temperature for the target pathogen and matrix to be tested prior to use. See individual assay instructions for further information.

6. DISPOSAL

Dispose of glass vials into an appropriate sharps waste container.



7. SAFETY

For use in laboratory facilities with qualified personnel trained for the handling of potentially pathogenic organisms. Training is recommended to first time users and can be provided by Solus Scientific Solutions Ltd. Using the method requires compliance with Good Laboratory Practices (refer to EN ISO 7218).

As a guide, the following precautions should be taken as a minimum:

- Protective clothing should be worn including lab coat, safety glasses, mask and gloves where appropriate.
- Do not pipette by mouth.
- Avoid contact with the skin.
- Do not inhale powder.
- Do not eat, drink or apply cosmetics in the laboratory.

Biological samples such as enrichments have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations on disposal of biological wastes. Wear appropriate protective equipment which includes but is not limited to: protective eyewear, face shield, clothing/laboratory coat, and gloves. All work should be conducted in properly equipped facilities utilizing the appropriate safety equipment (e.g. physical contaminant devices). Individuals should be trained in accordance with applicable regulatory and company/institution requirements before working with potentially infectious materials. All enrichment broths should be sterilized following any culture based confirmatory steps through heat denaturation by autoclaving at 121°C for 15 min.

8. PRECAUTIONS

- Optimum sensitivity and specificity will be reduced if contents are modified or not stored under the recommended conditions.
- Avoid microbial contamination of opened bottles.
- Do not use for diagnostic purposes with medical specimens.

9. MSDS INFORMATION

Material safety data sheets (MSDS) are available for this product on request.

Copyright Information

This document, including all photographs and illustrations, contains proprietary information that is protected by copyright. No part of this publication may be reproduced in any form whatsoever or translated into any language without the prior, written permission of PerkinElmer. ©2020 PerkinElmer, Inc. All rights reserved.

Trademarks

PerkinElmer is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

For more information visit www.solusscientific.com

PerkinElmer, Inc
940 Winter Street,
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs