



3050 Spruce Street
Saint Louis, Missouri 63103 USA
Telephone 800-325-5832 • (314) 771-5765
Fax (314) 286-7828
email: techserv@sial.com
sigma-aldrich.com

Product Information

Formalin Neutralizer Kit

Catalog Number **FN0010**
Store at Room Temperature

TECHNICAL BULLETIN

Product Description

The health hazards associated with formaldehyde, or formalin, are well known¹. Disposal of formalin through licensed waste facilities can be time consuming and expensive. Federal regulations² allow a user to treat their own waste without obtaining a permit and without being regulated as a Treatment, Storage, or Disposal facility. The Formalin Neutralizer Kit does not add organics or heavy metals to the neutralized product, making it waste drain disposable (state and local regulations prevail). The kit will neutralize 10 liters of 10% neutral buffered formalin.

Components

Formalin Neutralizer, Catalog Number F5429
Formalin Neutralizer Test Reagent, Catalog Number . F5679

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

The use of Personal Protective Equipment (PPE) and adequate ventilation are required when handling and mixing the formalin waste with the Formalin Neutralizer.

Storage/Stability

Keep tightly closed. The Formalin Neutralizer should be stored at Room Temperature. Upon receipt of the kit, it is recommended that the Formalin Neutralizer Test Reagent be stored at 2-8 °C.

Preparation Instructions

The Formalin Neutralizer is ready to use. Warm the Formalin Neutralizer Test Reagent to room temperature before use.

Procedure

Solution to be Treated	Quantity of Solution to be Treated	Formalin Neutralizer Required
1% formaldehyde	1 gallon	2 level scoops*
10% Neutral Buffered Formalin	1 liter	2 level scoops*
10% Neutral Buffered formalin	1 gallon	8 level scoops*

* Use scoop provided in kit

1. Add Formalin Neutralizer as indicated in the table above and mix until dissolved.
2. Wait at least 10 minutes.
3. Test neutralized formalin solution.
 - a. Add 100 µl of neutralized formalin solution to 100 µl of Formalin Neutralizer Test Reagent.
 - b. Cap the tube – invert 3 times to mix thoroughly
 - c. Wait 5 minutes for color to develop.
 - d. Test solutions which are colorless, light or medium violet or pink, have been neutralized.

Note: Colorless solutions have a formaldehyde content <0.001%. Solutions that yield a deep pink or burgundy color after treatment have a formaldehyde content >0.05% and should be retreated. Repeat Steps 1 to 3.

Run a positive control using untreated formalin waste solution – see Step 3 above. Solution should turn a deep pink to deep burgundy. If solution does not turn a deep burgundy, contact Sigma Technical Service for assistance.

References

1. United States Code of Federal Regulations. 29 CFR 1910.1048 (2003)
2. United States Code of Federal Regulations. 40 CFR 261.5 (2004)

RK,PHC 08/05-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.