

Product Information

Phenol Solution

P4557

Storage Temperature: 2–8 °C

Synonym: Equilibrated Phenol

Product Description

The product is a clear liquid capped with a layer of Tris buffer.

- pH of phenolic phase: 6.7 ± 0.2
- pH of phenolic phase: 7.9 ± 0.2 (after the addition of Equilibration Buffer)
- This preparation may be used in the extraction of nucleic acids.

Components

The product is supplied with Equilibration Buffer (B5658).

Precautions and Disclaimer

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

Preparation Instructions

For many applications, this product can be used without the addition of the Equilibration Buffer. For applications requiring a higher pH, such as the isolation of large intact genomic DNA, addition of the Equilibration Buffer (B5658) is recommended. Add the entire contents of the small bottle of Equilibration Buffer to the large bottle of Phenol solution. Mix gently and allow the phases to separate before use, 2–4 hours. This will adjust the pH of the Phenol Solution from $\text{pH } 6.7 \pm 0.2$ to 7.9 ± 0.2 .

Storage/Stability

Equilibrated Phenol is packaged under an inert gas and is for use if the phenolic phase is colorless.

Do not use if the phenolic phase has turned to a yellow or brown color. After the addition of Equilibration Buffer, the product may be stored for 6 months at 2–8 °C.

For extended storage of the product after opening, either with or without the addition of Equilibration Buffer, –20 °C with protection from light is recommended.

References

1. Sambrook, J., Fritsch, E. F., & Maniatis, T. (1989). *Molecular cloning: a laboratory manual* (No. Ed. 2, pp. xxxviii+-1546).

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page at SigmaAldrich.com/techservice.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/offices.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck and Sigma-Aldrich are trademarks of Merck or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2008-2026 Merck and/or its affiliates. All Rights Reserved.
P4557pis Rev 04/26

