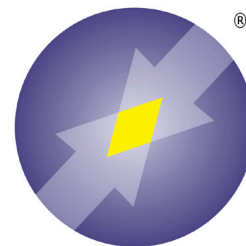


Kathon CG Preservative

Catalog Number X129-25ML



ARBOR
ASSAYS

FEATURES

- A broad spectrum preservative suitable for HRP containing materials
- Does **NOT** contain mercury or azide
- Stable solution at room temperature

INTRODUCTION

Kathon CG is an EPA-registered formaldehyde-free preservative for biological applications. This preservative displays exceptional antimicrobial activity against gram-positive and gram-negative bacteria, yeasts, and molds in a variety of solutions extending their shelf life and delaying decomposition.

It displays no inhibitory effect on peroxidase enzymes and, as such, is an ideal preservative for buffers and solutions that contain or will come in contact with HRP-labeled antibodies, conjugates and with peroxidase containing solutions. Typical concentrations for effective preservative action are $\leq 0.1\%$.

FORM:	Aqueous solution
VOLUME:	25 mL
STORAGE:	Room temperature
STABILITY:	9 months
USES:	As a preservative for aqueous buffers and solutions.
COUNTRY OF ORIGIN:	USA

FOR RESEARCH USE ONLY