

6033X-75**Microscopy****Harleco® Fuchsin, Carbol Solution Kinyoun**

AFB Stain



In Vitro Diagnostic Medical Device

Intended Use

Harleco® Fuchsin, Carbol Solution Kinyoun is used to stain cultures or specimens suspected to contain Mycobacteria (Acid-Fast Bacteria, or AFB). It serves the purpose of investigating sample material of human origin.

Principle

The Fuchsin, Carbol Solution Kinyoun is the conventional "cold staining" method. Positive staining reactions are evidence of Mycobacteria in the specimen; however, further analysis by culturing methods are required to confirm identification. All specimens submitted for determination of possible mycobacterial infection should be examined for the presence of acid-fast *bacilli*. A major characteristic of mycobacteria is that they are "acid-fast" and once stained with an aniline dye, basic fuchsin, they are difficult to decolorize thus retaining the red color even when exposed to an acid alcohol rinse.

Sample material

Staining may be performed on smears prepared from specimens or samples containing mixed flora or pure cultures. It is recommended that smears be placed low on the slide to maximize reagent use. Specimens may be smeared directly on clean unused slides. Smears are allowed to air dry and then passed through a low flame two to three times to heat fix.

Caution: DO NOT OVER FIX WITH EXCESSIVE HEAT!

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the US and Canada.

Reagent

Cat. No. 6033X 1 L
Harleco® Fuchsin, Carbol Solution Kinyoun

Also required:

Cat. No. A3179 Acid Alcohol, 70% 1 L
Cat. No. 1808-50ML Methylene Blue Solution 50 mL

Sample preparation

The sampling must be performed by qualified personnel. All samples must be clearly labeled. Suitable instruments must be used for collecting and preparing samples. Follow the manufacturer's instructions for application/use.

Manual Staining Procedure

1. Using slide holder or individual slides, submerge slide in the Carbol Fuchsin Solution for 5 minutes.
2. Rinse gently with deionized water.
3. Remove most of the stain by using the Acid Alcohol (Decolorizer) Solution until the washings are clear.
4. Rinse gently with tap water.
5. Immerse the slide in Methylene Blue Solution for 10-20 seconds.
6. Rinse gently with tap water.
7. Allow to air dry.

Results

Acid-fast mycobacteria	Red
Non-acid fast mycobacteria	Blue
Background	Blue

Diagnostics

Diagnoses are to be made only by authorized and trained personnel. Valid nomenclature must be used. Further tests must be selected and implemented according to recognized methods. Suitable controls should be conducted with each application.

Storage

15-30 °C

Shelf-life

The Fuchsin, Carbol Solution Kinyoun for microscopy can be used until the stated expiry on the packaging.

After first opening of the bottle, the contents can be used up to the stated expiry date when stored at 15-30 °C.

The bottles must always be kept tightly closed.

Additional instructions

For professional use only.

The application must be carried out by qualified personnel only.

National guidelines for work safety and quality assurance must be followed.

Microscopes equipped according to the standards must be used.

Protection against infection

Effective measures must be taken to protect against infection in line with laboratory guidelines.

Instructions for disposal

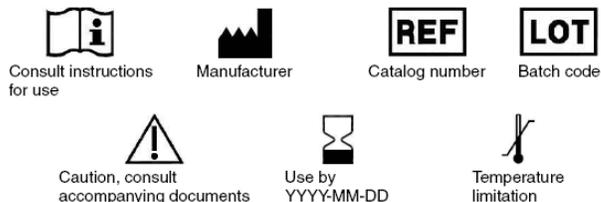
The package must be disposed of in accordance with the current disposal guidelines. Used solutions and solutions that are past their shelf-life must be disposed of as special waste in accordance with local guidelines.

Auxiliary reagents

Cat. No. 64969	Harleco® Krystalon™ Mounting Medium	50 mL, 500 mL
Cat. No. 104699	Immersion oil for microscopy	100 mL, 500 mL
Cat. No. 103699	Immersion oil acc. To ISO 8036, halogen-free	100 mL

Hazard classification

Cat. No. 6033X Please observe the hazard classification printed on the label and the information given in the safety data sheet. The safety data sheet is available on the website and on request.



Literature

1. Conn's Biological Stains: A Handbook of Dyes, Stains and Fluorochromes for Use in Biology and Medicine, 10th Edition, (ed. Horobin, R.W., and Kiernan, J.A.) Bios, 2002

Harleco® is a registered trademark of Merck KGaA, Darmstadt, Germany.

Krystalon™ is a trademark of Merck KGaA, Darmstadt, Germany

Status: 2020-07-10

20499381



EMD Millipore Corporation
400 Summit Drive
Burlington, MA 01821, USA
Tel. +1-978-715-4321