

Thermo Scientific Xylene Instructions for Use

For in vitro diagnostic use.

For the preparation of pathology specimens.

Thermo Scientific™ Xylene is specifically designed for tissue processing and staining of histological and cytological specimens. The product is a clear, colorless reagent that is highly quality controlled. Gas chromatography techniques are performed on each batch of Xylene to ensure chemical consistency and to identify any impurity that may affect tissue processing and/or staining. Benzene, a known carcinogen, is a contaminant of all xylene (xylol). Our quality assurance program ensures the lowest level of benzene. Precautions should be taken when handling Xylene. Protective gloves should be worn and work should be performed in a ventilated area. Xylene can be used with both open and closed tissue processors. It is also compatible with manual staining procedures and automatic stainers. Xylene is a flammable product and must be stored in a flammable fire cabinet.

Tissue Processing Instructions For Use

Xylene has shown best results on closed tissue processors when heat is not used during the normal clearing process. Vacuum has been viewed as an asset during all phases of tissue processing, including clearing. A typical tissue processing schedule would include two stations and the time in each Xylene station should be 1 hour for a multiple tissue type/thickness processing run. Xylene is also recommended for use in the cleaning cycle of a closed tissue processor.

The laboratory should develop a product rotation and replacement schedule that adheres to the policies of their department.

Protocols

The following overnight tissue processing schedule is provided as an example for reference purposes.

Station	Solution	Time
1	10% Neutral Buffered Formalin	holding
2	10% Neutral Buffered Formalin	1 hour
3	Pen-Fix™ or 80% Alcohol	40 minutes
4	95% Alcohol	40 minutes
5	95% Alcohol	40 minutes
6	100% Alcohol	40 minutes
7	100% Alcohol	40 minutes
8	100% Alcohol	40 minutes
9	Xylene	1 hour
10	Xylene	1 hour
11	Paraffin	1 hour
12	Paraffin	1 hour

Note: This procedure may not fit every situation. Modifications may be necessary.

The following biopsy (less than 2 mm in thickness) tissue processing schedule is provided as an example for reference purposes. Tissues are assumed to be fixed. If not, stations 1 and 2 should utilize 10% Neutral Buffered Formalin for a minimum of 30 minutes each.

Station	Solution	Time
1	10% Neutral Buffered Formalin	(30 minutes)
2	10% Neutral Buffered Formalin	(30 minutes)
3	Pen-Fix or 80% Alcohol	10 minutes
4	95% Alcohol	10 minutes
5	95% Alcohol	10 minutes
6	100% Alcohol	10 minutes
7	100% Alcohol	10 minutes
8	100% Alcohol	10 minutes
9	Xylene	15 minutes
10	Xylene	15 minutes
11	Paraffin	20 minutes
12	Paraffin	20 minutes

Note: This procedure may not fit every situation. Modifications may be necessary.

Staining Instructions For Use

It is recommended that three stations of Xylene for 3 minutes each be used for deparaffinization to ensure all of the paraffin is removed from the tissue section.

After staining and dehydration is completed, slides should be cleared in three stations of Xylene for 1 minute each before coverslipping. This will ensure complete clearing, which results in maximum slide clarity, and readies the slide for coverslipping.

The laboratory should develop a product rotation and replacement schedule that adheres to the policies of their department.

Protocol

The following H&E staining schedule is provided as an example for reference purposes.

Station	Solution	Time
1	Xylene	3 minutes
2	Xylene	3 minutes
3	Xylene	3 minutes
4	100% Alcohol	1 minute
5	100% Alcohol	1 minute
6	100% Alcohol	1 minute
7	95% Alcohol	1 minute
8	Rinse in running tap water	Briefly
9	Deionized or distilled water	Rinse
10	Hematoxylin	Chosen Time
11	Running tap water	Rinse off excess stain
12	Acid Rinse	Chosen Time
13	Rinse in running tap water	30 seconds (agitate)
14	Bluing Reagent	1 minute
15	Rinse in running tap water	1 minute
16	95% Alcohol	Rinse
17	Eosin-Y	Chosen Time
18	100% Alcohol	1 minute
19	100% Alcohol	1 minute
20	100% Alcohol	1 minute
21	Xylene	1 minute
22	Xylene	1 minute
23	Xylene	1 minute

Note: This procedure may not fit every situation. Modifications may be necessary.

Warnings and Precautions

See Safety Data Sheets for warnings and precautions, as well as EUH code definitions.

See container label for warnings and precautions.

Order Information

Product	Size	Qty.	REF
Richard-Allan Scientific™ Xylene	1 gal. (3.79 L) bottle	4/cs.	6601
Richard-Allan Scientific™ Xylene	1.32 gal. (5 L) bottle	4/cs.	6615
Syntri Safeguard™*			
Richard-Allan Scientific™ Xylene	1.32 gal. (5 L) bottle	4/cs.	6615SS
Richard-Allan Scientific™ Xylene	55 gal. (208 L) drum	Ea.	6655
Richard-Allan Scientific™ Xylene (ACS)	1 gal. (3.79 L) bottle	4/cs.	9900-1
Richard-Allan Scientific™ Xylene (ACS)	5 gal. (18.9 L) metal pail	Ea.	9900-5
Richard-Allan Scientific™ Xylene (ACS)	55 gal. (208 L) drum	Ea.	9900-55
Shandon™ Xylene	1.06 gal. (4 L) bottle	Ea.	9990501
Shandon™ Xylene	2.64 gal. (10 L) bottle	2/cs.	9990502
Shandon™ Xylene	55 gal. (208 L) drum	Ea.	9990503

*For further instructions on how to use the Syntri Safeguard products please consult the Thermo Scientific Revos user's manual

Anatomical Pathology



IS8126 R10/18



Richard-Allan Scientific
Subsidiary of Thermo Fisher Scientific
4481 Campus Drive, Kalamazoo, MI 49008
Phone: 1 (800) 522-7270
Phone: + 1 (269) 544-5600
www.thermofisher.com/pathology



Thermo Shandon Limited
Subsidiary of Thermo Fisher Scientific
Tudor Road, Manor Park,
Runcorn, Cheshire WA7 1TA, UK
Phone: +44 (0) 1928 534 050
Fax: + 44 (0) 1928 534 049



ThermoFisher
SCIENTIFIC
The world leader in serving science