

Thermo Scientific Richard-Allan Scientific 10% Neutral Buffered Formalin Instructions for Use

For in vitro diagnostic use.

For fixation of pathology specimens.

Thermo Scientific™ Richard-Allan Scientific™ 10% Neutral Buffered Formalin is designed to be used at full strength and should not require any additives. It is used as a primary fixative.

Specimen Preparation

1. Tissue specimens should be placed in 10% Neutral Buffered Formalin immediately after removal from the body. Prompt fixation will prevent autolysis, putrefaction, and other undesirable cellular changes.
2. When grossing, tissue specimens should be no thicker than 5mm to facilitate optimal fixation.
3. Containers should be large enough to accommodate the specimen and to allow a ratio of 15 to 20 times the volume of 10% Neutral Buffered Formalin to the volume of the specimen.

Procedure

1. 10% Neutral Buffered Formalin is commonly used in the grossing room as a holding solution until the tissue specimens are ready to be placed onto the tissue processor. This solution should be replaced daily.
Warning: Use this product in a well ventilated area and use protective gloves and other precautions when handling 10% Neutral Buffered Formalin.
2. Use 10% Neutral Buffered Formalin in the first two stations of the tissue processor to ensure adequate fixation of tissue sections. One hour immersion time in each station is generally used in conventional tissue processing. 10% Neutral Buffered Formalin can be used on all tissue processors.
3. Tissue specimens can be stored in containers filled with 10% Neutral Buffered Formalin indefinitely. Specimens will not over-harden or be prone to excessive shrinkage from prolonged storage.
4. If used in enclosed tissue processors, it is advisable to follow the manufacturer's instructions regarding flushing and housekeeping procedures.

For 2.5 and 5 Gallon Cube

The 2.5 and 5 gallon cubes with spigot are used to easily dispense 10% Neutral Buffered Formalin into tissue processor reagent vessels or specimen containers. Simply remove the cap and replace with the supplied spigot and fasten tightly.

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
10% Neutral Buffered Formalin	1 gal. (3.79 L) bottle	4/cs.	5701
10% Neutral Buffered Formalin	5 gal. (18.9 L) cube	Ea.	5705
10% Neutral Buffered Formalin	5 gal. (18.9 L) cube	Ea.	5705C
10% Neutral Buffered Formalin	1.32 gal. (5 L) bottle	4/cs.	5735
Syntri Safeguard™*			
10% Neutral Buffered Formalin	1.32 gal. (5 L) bottle	4/cs.	5735SS
Syntri Safeguard™*			
10% Neutral Buffered Formalin	1.32 gal. (5 L) bottle	4/cs.	5735SSC
10% Neutral Buffered Formalin	2.5 gal. (9.48 L) cube	Ea.	5725
10% Neutral Buffered Formalin	2.5 L	4/cs.	5745
10% Neutral Buffered Formalin	2.5 L	4/cs.	5745C
10% Neutral Buffered Formalin	55 gal. (208 L) drum	Ea.	5755
High Purity 10% Neutral Buffered Formalin	1 gal. (3.79 L) bottle	4/cs.	9400-1
High Purity 10% Neutral Buffered Formalin	5 gal. (18.9 L) cube	Ea.	9400-5
High Purity 10% Neutral Buffered Formalin	55 gal. (208 L) drum	Ea.	9400-55

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

