

Thermo Scientific Richard-Allan Scientific Acetone Instructions for Use

For in vitro diagnostic use.

For the preparation of pathology specimens.

Thermo Scientific™ Richard-Allan Scientific™ Acetone is specifically designed for laboratory procedures. The product is a clear, colorless reagent that is highly quality controlled. Acetone is a known toxin. Ingestion may result in blindness or death. Precaution should be taken when handling Acetone. Protective gloves should be worn and work should be performed in a ventilated area. Acetone is a highly flammable product and must be stored in a flammable fire cabinet.

Instructions For Use

Fixation

Acetone is a non-additive, protein coagulating fixative that is used when the demonstration of enzymes, (e.g. acid and alkaline phosphatase) is required. Fixation with Acetone is done rapidly in the refrigerator and most of the dehydration is accomplished at the same time. It is also used for fixation of brain tissue when staining techniques for rabies diagnosis is required. Acetone is frequently used on frozen sections of tissue to be stained by immunohistochemical techniques. Acetone is a very rapid acting fixative that causes extreme shrinkage, distortion and over-hardening; it is recommended for preservation of special tissue components.

Dehydration

Acetone is a very rapid acting dehydrant and causes excessive shrinkage of the tissue. Acetone is very volatile therefore maintaining proper solution levels on an open tissue processor is difficult. It is also very flammable and should not be used on a closed tissue processor.

Clearing

Acetone has a very low boiling point and will boil off when placed into paraffin at 58° C or higher. Acetone is also very effective in clearing and identifying lymph nodes in mesentery tissue.

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
Acetone	1 gal. (3.79 L)	4/cs.	9011
Acetone	5 gal. (18.9 L) metal pail	Ea.	9005
Acetone (ACS)	1 gal. (3.79 L) bottle	4/cs.	9000-1
Acetone (ACS)	5 gal. (18.9 L) metal pail	Ea.	9000-5
Acetone (ACS)	55 gal. 208 L drum	Ea.	9000-55

