

**Thermo Scientific Richard-Allan Scientific Mounting Media, Non-Aqueous
Instructions for Use**

For in vitro diagnostic use.

For use as a permanent mounting medium in slide preparation.

Mounting media are aromatic based adhesives for use with glass coverslips. They are clear and colorless and will not discolor with age. The formulas contain an antioxidant to inhibit stain fading. Their refractive index is 1.495 +/- 0.005 and they are soluble in xylene, toluene, and some aliphatic hydrocarbon based clearing reagents.

Thermo Scientific™ Richard-Allan Scientific™ Mounting Medium is a toluene-based rapid drying mounting medium. It is miscible with all clearing agents including xylene substitutes. This mounting medium is optically clear and will not cause fading of stains even after prolonged storage.

Thermo Scientific™ ClearVue™ Mountant is a toluene-based rapid drying mounting medium. It is miscible with all clearing agents including xylene substitutes. This mounting medium is optically clear and will not cause fading of stains even after prolonged storage. ClearVue mountant is packaged specifically for use with the Thermo Scientific™ ClearVue™ coverslipper.

Thermo Scientific™ ClearVue™ Mountant XYL is a xylene-based rapid drying mounting medium. It is miscible with all clearing agents including xylene substitutes. This mounting medium is optically clear and will not cause fading of stains even after prolonged storage. ClearVue mountant is packaged specifically for use with the Thermo Scientific™ ClearVue™ coverslipper.

Thermo Scientific™ Richard-Allan Scientific™ Cytoseal™ 60 is a toluene-based mounting medium formulated from acrylic resins and will not crack or discolor with age. The addition of antioxidants inhibits fading or yellowing of stained specimens. This is a low-viscosity medium that dries quickly and allows for rapid, even spreading which virtually eliminates air bubbles.

Thermo Scientific™ Richard-Allan Scientific™ Cytoseal™ 280 is a toluene-based mounting medium formulated from acrylic resins and will not crack or discolor with age. The addition of antioxidants inhibits fading or yellowing of stained specimens. This is a high-viscosity medium for applications where minimal spreading is desired or without a cover glass.

Thermo Scientific™ Richard-Allan Scientific™ Cytoseal™ XYL is a xylene-based mounting medium that contains an antioxidant to inhibit the fading or yellowing of stained specimens. It is a rapid drying medium allowing microscopic examination of slides soon after application.

Instructions For Use

1. Lay a coverslip on an absorbent surface.
2. Place a drop or two of mounting medium toward the edge of the coverslip.
3. Remove a stained slide from clearing reagent.
4. Quickly wipe the back of the slide.
5. Invert the slide so that the specimen side faces down.
6. Lower the slide until it reaches the coverslip, starting at one long edge. Gently press until the mounting medium spreads over the entire surface to be joined.
7. Turn over the slide/coverslip assembly.
8. If necessary, drain extra medium by standing the slide on its edge on an absorbent towel.
9. Gently press out any noticeable air bubbles with a pair of forceps.

Note: For Thermo Scientific Richard-Allan Scientific Mounting Medium catalog #4111, 4112, 4211 and 4212, do not use with picric acid counterstains; the stained slide will fade.

Removing a Coverslip

It is sometimes necessary to remove a coverslip. This is usually done so that a slide can be de-stained and re-stained. In this way a slide that was stained improperly can be corrected, or the original stain can be removed to make room for the application of a special stain.

To remove a coverslip do one of the following:

1. Soak the slide in xylene until the coverslip falls off.
 2. Place the slide in a 60° C oven for 3 to 4 hours.
 3. Place the slide in a freezer for several minutes.
- Do not try to force a coverslip off a slide. Damage to the tissue section will occur.

Automated Coverslipping

Follow directions for use with your automated coverslipping instrument.

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
Mounting Medium	4oz. (118 mL)	6/cs.	4112
Mounting Medium	16oz. (473 mL)	2/cs.	4111
ClearVue Mountant	16oz. (473 mL)	2/cs.	4211
ClearVue Mountant XYL	16oz. (473 mL)	2/cs.	4212
Cytoseal 60	4oz. (118 mL)	6/cs.	8310-4
Cytoseal 60	16oz. (473 mL)	6/cs.	8310-16
Cytoseal 280	4oz. (118 mL)	6/cs.	8311-4
Cytoseal XYL	4oz. (118 mL)	6/cs.	8312-4