

## Thermo Scientific DPX Mounting Media Instructions for Use

**For in vitro diagnostic use.  
For use as a permanent, resin-based mounting medium  
in histological and cytological preparations.**

### Intended Use

The Thermo Scientific™ DPX Mounting Media is used as a synthetic mounting medium in histology and cytology.

### Warnings and Precautions

#### For in vitro diagnostic use

Reasonable care should be taken when using all laboratory reagents. Use with adequate ventilation.

Contains xylene 60 - 70%.

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

### Storage

Keep in a cool, dry well ventilated area. Keep containers tightly closed. Store in correctly labelled containers. Keep away from direct sunlight.

### Principles of the Method

DPX is suitable for all staining techniques which are compatible with the use of alcohol and aromatic (xylene/ toluene) clearing agent. DPX is clear and colorless and will not discolor with age. The formula contains an anti-oxidant to inhibit stain fading, and is completely soluble in xylene and toluene. DPX is dispensed onto a stained specimen slide in liquid form. The evaporation of the carrier solvent results in the hardening of the DPX to form a solid clear film. As DPX has a reflective index similar to glass, the specimen can be viewed under a microscope.

### Instructions for Use

1. Settle 8 hours prior to use, to allow any micro bubbles formed during transport to settle out.
2. Lay a coverslip on an absorbent surface.
3. Manually, using a glass rod or pipette, drop approximately 1 drop (0.5 mL) of DPX toward the edge of the coverslip.
4. Remove a stained slide from the clearing agent.  
**Note:** The stained slide should be well drained of clearing agent prior to mounting.
5. Invert the stained slide so that the slide faces down.
6. Lower the stained slide until it reaches the coverslip.
7. Starting at one long edge, gently press the stained slide until the DPX spreads over the surface area of the coverslip.
8. Turn over the stained slide and coverslip assembly and gently press out any noticeable air bubbles with a pair of forceps. Any excess DPX should be removed using an absorbant towel.
9. Allow the stained slide and coverslip assembly to dry in a horizontal position for one hour, prior to moving or viewing under a microscope to avoid movement of the coverslip.  
**Note:** The use of excessive heat to dry the slides is not recommended.
10. The coverslip and DPX can be removed by soaking the mounted slide in a corresponding clearing agent until the coverslip and mounting medium can be removed.
11. Allow the mounted slides to dry for 48 hours prior to filing. This ensures that slides are thoroughly dry and prevents slides sticking together.
12. Stained slides should be stored in a dark environment, away from sources of bright light, and checked periodically to confirm the DPX remains clear; with no sign of the DPX drying back or peeling off the slide.
13. DPX is a ready to use product, any modification is done so at the user's own risk.

#### Follow directions for use with your automated coverslipping instrument.

**Note:** Thermo Scientific has not validated the protocol illustrated here and takes no responsibility for its use. Customer should always validate protocols before placing any reliance on them.

### Order Information

Product	Size	REF
DPX	1 L	LAMB/DPX

