

# Compatibility of Mounting Media on Thermo Scientific Nunc Permanox Slides

We use two types of surfaces for Thermo Scientific Nunc Lab-Tek Chamber Slide System, glass and plastic. One of the plastics used is Thermo Scientific Nunc Permanox, which is a special, oxygen permeable polymer.

This polymer has excellent optical qualities and is resistant to many commonly used solvents such as acids, alkalis, organic solvents such as ethanol and some

hydrocarbons such as ethylene glycol.

However, Nunc™ Permanox™ is not chemically compatible with any organic reagent containing solvents such as xylene or toluene. These solvents will warp the slides.

The following table lists some aqueous mounting media that can be used on the Lab-Tek™ Permanox slides for various applications. Permanox slides were

coverslipped with the different mounting media listed in the table. Mounting medium/slide compatibility was determined by the degree of warpage after one week at room temperature. The mounting medium application is as recommended by manufacturer's product information. Stain/mounting medium compatibility must be optimized for each application.

Media Company Product Name	Product Description	Media Application	Compatible with Permanox Plastic
<b>Thermo Scientific Shandon Permafluor</b>	<b>Aqueous</b> <ul style="list-style-type: none"> <li>• Permanent</li> <li>• Recommended storage: 2-8°C</li> <li>• Use room temperature</li> </ul>	<b>Fluorescence microscopy</b> <ul style="list-style-type: none"> <li>• Enhances FITC intensity and reduces fading</li> </ul>	Yes
<b>Thermo Scientific Shandon Immu-Mount</b>	<b>Aqueous</b> <ul style="list-style-type: none"> <li>• Permanent</li> <li>• Recommended storage: room temperature</li> </ul>	<b>Immunocytochemical staining with frozen sections</b>	Yes
<b>Biomedica Gel Mount</b>	<b>Aqueous</b> <ul style="list-style-type: none"> <li>• Permanent</li> <li>• Recommended storage: room temperature</li> </ul>	<b>Immunofluorescence</b> <ul style="list-style-type: none"> <li>• Use in conjunction with fluorescein, rhodamine, Texas Red and Phycobiliproteins</li> </ul>	Yes
<b>Biomedica Crystal Mount</b>	<b>Aqueous</b> <ul style="list-style-type: none"> <li>• Permanent</li> <li>• Recommended storage: room temperature</li> <li>• No coverslips required</li> <li>• Heat slide in oven 70-80°C, 10 minute</li> </ul>	<b>Immunocytochemical staining (AEC, BCIP, NBT, DAB)</b> <ul style="list-style-type: none"> <li>• Immunofluorescence</li> <li>• Compatible with common fluorochromes</li> </ul>	No* *Slides can warp if heated
<b>Vector Laboratories Vectashield</b>	<b>Aqueous</b> <ul style="list-style-type: none"> <li>• Not Permanent</li> <li>• Recommended storage: 4°C in dark</li> </ul>	<b>Immunofluorescence</b>	Yes

[www.thermoscientific.com](http://www.thermoscientific.com)

© 2010 Thermo Fisher Scientific Inc. All rights reserved. "Gel Mount" and "Crystal Mount" are registered trademarks of Biomedica; and "Vectashield" is a registered trademark of Vector Laboratories. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

**Austria**

+43 1 801 40 0

**Belgium**

+32 53 73 42 41

**China**

+86 21 68654588

**Denmark**

+45 4631 2000

**France**

+33 2 2803 2180

**Germany**

+49 6184 90 6940

**India**

+91 22 6716 2200

**Italy**

+39 02 02 95059 or  
434-254-375

**Japan**

+81 3 3816 3355

**Netherlands**

+31 76 571 4440

**Nordic/Baltic  
countries**

+358 9 329 100

**North America**

+1 585-586-8800

**Russia/CIS**

+7 (812) 703 42 15

**Spain/Portugal**

+34 93 223 09 18

**South America**

+1 585 899 7298

**Switzerland**

+41 44 454 12 12

**UK/Ireland**

+44 870 609 9203

**Other Asian countries**

+852 2885 4613

**Countries not listed**

+49 6184 90 6940 or  
+33 2 2803 2180

TILSPNUNCTN15 0610