

# VectaMount<sup>®</sup> Mounting Medium

<b>Cat. No.</b>	H-5000
<b>Storage</b>	Room Temperature
<b>Description</b>	Optically clear and odorless formula for permanently preserving histochemical stains or precipitable enzyme substrates in tissue sections or cell preparations.

VectaMount Mounting Medium is a non-flammable, non-hazardous, and odorless formula for permanently preserving precipitable enzyme substrates in tissue sections or cell preparations. VectaMount Mounting Medium retains the color and intensity of preparations attained with commonly used enzyme substrates. It is pale yellow in color, optically clear, and contains no toluene or xylene. VectaMount Mounting Medium dries clear with an ideal refractive index.

VectaMount Mounting Medium is compatible with most clearing agents\*, including aliphatic hydrocarbon- (e.g. d-limonene) or aromatic hydrocarbon- (toluene or xylene) based clearing agents.

## Kit Components

Product Name	Volume
VectaMount Mounting Medium	60 mL

\* For most xylene-free clearing agents such as Clear-Rite 3, incubation times of at least 3 x 3-5 minutes is recommended.

## Guidelines for Use

Following dehydration and clearing of the tissue section or cell preparation on a slide, dispense a small amount of VectaMount Mounting Medium on or adjacent to the section using a pipette (plastic or glass) or glass rod. As the coverslip is applied, the medium will disperse over the entire coverslipped area.

Allow the preparation to dry. This may take one to several hours, depending on the temperature, humidity, type of clearing agent used, and amount of mounting medium applied. Drying can be accelerated by incubating the slide in a 60-75 °C oven for 15-30 minutes. Allow the slide to cool before storage.

If necessary, the coverslip can be removed by incubating the slide in xylene or a d-limonene-based clearing agent.

Keep container tightly sealed when not in use. Some cloudiness of the solution may be present but will have no adverse affect on the appearance of the mounted section.

Detailed product listings, specifications, protocols and additional information is available on our website: [vectorlabs.com](http://vectorlabs.com)