

## VectaMount® AQ Mounting Medium

**Cat. No.:** H-5501

**Storage:** Room Temperature

A hardsetting mounting medium developed for use with enzymatic substrates, whose reaction products are soluble in alcohol or other organic solvents

### Principle:

VectaMount AQ Mounting Medium is a hard-setting aqueous-based mounting medium developed for use with enzymatic substrates, whose reaction products are soluble in alcohol or other organic solvents. VectaMount AQ Mounting Medium is also compatible with most peroxidase and alkaline phosphatase enzyme substrates.

### COMPONENTS

<u>Product Name</u>	<u>Volume</u>
VectaMount AQ Mounting Medium	60 ml



### Guidelines for Use:

To avoid bubbles, do not shake bottle. Invert bottle and dispense drops, keeping constant pressure on the bottle. If there are any bubbles in the bottle tip, dispense and discard several drops before use.

Squeeze one or two drops of VectaMount AQ Mounting Medium onto the stained preparation. Coverslip and allow mounting medium to disperse over the entire section. Recap bottle when not in use.

VectaMount AQ Mounting Medium dries completely after slides are coverslipped. Sections may be viewed immediately, but generally 10-15 minutes are required for the coverslip to become immobilized and several hours for VectaMount AQ Mounting Medium to completely harden.

If air bubbles are present, or if a drying artifact develops on the slide during prolonged storage, coverslips can be removed and the sections remounted. Mounted coverslips can be easily removed by soaking in buffer (PBS or TTBS). Mounted slides can be stored at room temperature for at least 1 year without the substrate fading.

**Note:** For immunofluorescent sections or cells, VECTASHIELD® Vibrance™ Antifade Mounting Media are recommended to prevent photobleaching.

*\* VectaMount AQ Mounting Medium may also be stored at 2-8 °C but warming to room temperature is recommended before use.*

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