



CERTIFICATE OF ANALYSIS

Product *VectaMount*[™] Mounting Medium
Catalog H-5501
Amount 60 ml
Lot No. W0908
Storage Conditions Room Temperature*

VectaMount[™] AQ Mounting Medium is a hard-setting aqueous-based mounting medium developed for use with enzymatic substrates such as AEC, whose reaction products are soluble in alcohol or other organic solvents. *VectaMount*[™] AQ is also compatible with most peroxidase and alkaline phosphatase enzyme substrates, such as DAB, BCIP/NBT, Vector[®] SG, Vector[®] Red, Vector[®] Blue.

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 onto the stained preparation. Coverslip and allow mounting medium to disperse over the entire section. Recap bottle when not in use.

VectaMount[™] AQ 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 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 TBS). Mounted slides can be stored at room temperature for at least 1 year without the substrate fading.

Note:

For immunofluorescent sections or cells, VECTASHIELD[®] Hard+Set[™] (H-1400, H-1500) products 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.