

VECTABOND[®] Reagent

Tissue Section Adhesion

Cat. No. SP-1800

Storage Room Temperature

Description The VECTABOND Reagent for Tissue Section Adhesion significantly increases the adherence of both frozen and paraffin-embedded tissue sections and cell preparations to glass slides by chemically modifying the surface of the glass with positive charges to form a highly adherent surface.

Components

Product Name	Volume
VECTABOND Reagent	7 ml

Guidelines for Use

For best results, we recommend that the entire contents be used at once. If a partially used bottle is to be kept, store the sealed bottle at room temperature and protect from moisture.

Use acetone-compatible containers such as glass staining dishes for these reagents.

Slide Preparation & Treatment

1. Place clean glass slides in metal slide racks.
2. Immerse slides in acetone for 5 minutes.
3. Immediately before use, prepare VECTABOND Reagent treatment solution by adding the entire contents (7 ml) of the bottle to 350 ml of acetone and stir well.
4. Remove slides from acetone, tapping rack several times to drain and immediately place in VECTABOND Reagent treatment solution for 5 minutes.
5. Remove rack from VECTABOND Reagent and drain excess solution from rack. Gently dip slide rack in deionized or distilled water 3-4 times to remove excess reagent (do not create bubbles), change the water after every 5 racks.
6. Remove rack of slides from water and drain excess water. Air dry slides thoroughly at room temperature or at 37 °C.

Detailed product listings, specifications, protocols and additional information is available on our website: vectorlabs.com