



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NB900-66726](http://www.novusbio.com/NB900-66726)

## NB900-66726 Protocol

### Immunohistochemistry-Paraffin protocol for Fluoromount (NB900-66726)

---

Mounting Procedure (NB900-66726): [https://www.novusbio.com/products/fluoromount\\_nb900-66726](https://www.novusbio.com/products/fluoromount_nb900-66726)

1. Remove the red tip from the bottle and cut the tip off with a sharp razor or scissors. This will help achieve smooth flow of Fluoromount and prevent formation of tiny bubbles.
2. Place the bottle upside down in a container before using. This will help clear Fluoromount of tiny bubbles.
3. Blot excess water from the slide without letting tissue dry. Make sure the tissue is wet prior to mounting.
4. Apply 2-3 drops of Fluoromount on the tissue section and gently apply a cover slip. Allow to dry for 30-45 minutes at room temperature. Seal edges of the coverslip with nail polish for long term storage.