



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NB900-66726

NB900-66726 Protocol

Removal of Fluoromount (NB900-66726)

Fluoromount is an aqueous-based mounting medium and can easily be removed by soaking slides in deionized water. Place slide in a beaker full of deionized water on a magnetic stirrer. Leave the slide stirring in water for few hours or overnight for complete removal of dry Fluoromount.