



Mounting Medium
Catalog Number: CTS011

Reagent and Storage Specifications:

Component	Permanent Aqueous Mounting Medium is ready to use as supplied (15 mL).
Storage	Store at room temperature.

Protocol:

1. Complete the staining and counterstaining.
2. Apply the Mounting Medium.
3. Coverslip slides.
4. Let dry at room temperature.