

## Preparing the Ultramark for Shipment

You must remove the **Filter Magazine** and attach the safety **Lock Plates** before shipping the Ultramark. **Failure to do so will result in serious damage to the instrument.**

