

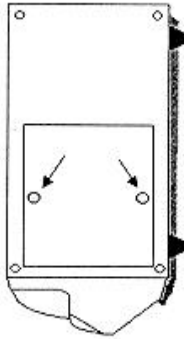
Important Information!

The Ultramark optical block has been locked for transportation and the Filter Magazine is not installed. Do not turn the instrument on before releasing the safety lock and installing the Filter Magazine.

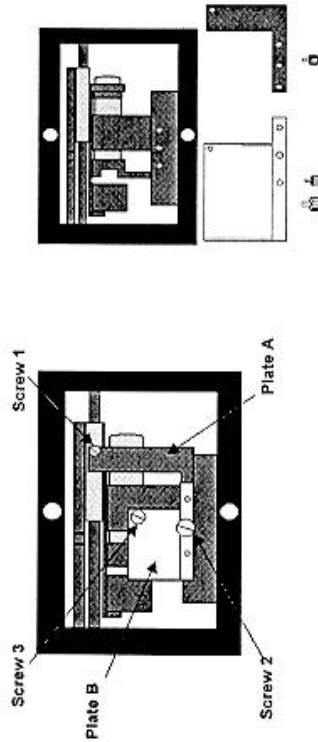
Follow the procedure below to unlock the instrument and install the Filter Magazine.

Releasing the Safety Lock

Remove the two thumb screws on the right side of the Ultramark and remove the side cover.

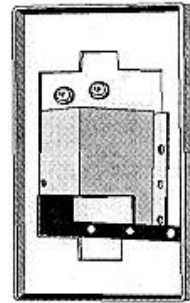


Inside the side panel, remove Screw 3 and Screw 2 to remove Lock Plate B. Remove Screw 1 to remove Lock Plate A.



NOTE: Screw 2 is used to install the Filter Magazine.

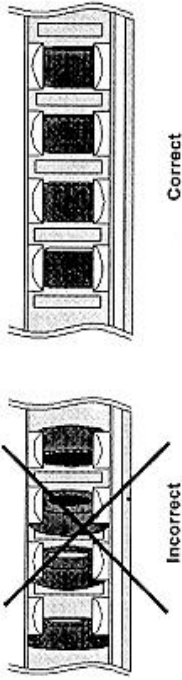
Keep the removed Lock Plates and Screws 1 and 3 in case you need to transport the instrument again. You can store them inside the box that contains the Filter Magazine.



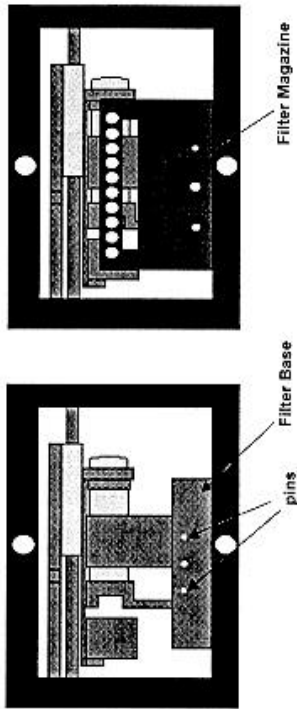
Installing the Filter Magazine

Remove the Filter Magazine from its packaging.

NOTE: Make sure that all the filters sit securely and properly in the Magazine clips before installation. Also, make sure that each filter in the installed Magazine is oriented with the numbered side facing the rear of the Ultramark and the shiny side facing the front.

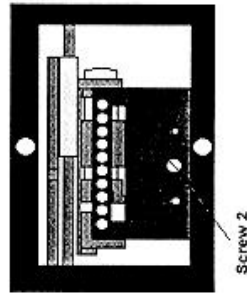


Insert the Filter Magazine into the base so that the pins in the base fit in the holes in the magazine.



Insert Screw 2 into the center hole in the Filter Magazine and tighten securely with a coin or screwdriver.

NOTE: The Filter Magazine must be securely fastened to the base to avoid serious problems.



Replace the side cover and fasten it with the thumb screws. The Ultramark is now ready to use.