

# Maintenance Checklist — Autopure LS

<b>Month:</b>																															
	Day																														
<b>Daily maintenance</b>	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31
Open the reagent cabinet and check the level of each reagent bottle, replace if necessary.																															
Visually check the reagent lines and ensure there are no air bubbles in any of the reagent lines.																															
Check for condensation in the centrifuge chamber and buckets. Wipe dry, if necessary.																															
Wipe any spills from the instrument worktable using a paper towel sprayed with cleaner.																															
Check the level of the waste bottle. Ensure the bottle has adequate room for additional Autopure LS runs.																															
<b>Optional:</b> Perform a Pump Dispense Test and verify that the correct amount of reagent was dispensed to all 8 tubes as described in the <i>Autopure LS User Manual</i> .																															
<b>Weekly maintenance</b>	Week 1		Week 2		Week 3		Week 4		Week 5																						
Wipe out the base of the waste tray using a paper towel sprayed with cleaner.																															
Shut down Windows. Turn off Autopure LS instrument. (Switch is on the left backside of the instrument.) Wait a few minutes. Turn the instrument on again.																															
<b>Monthly maintenance</b>	Date performed																														
Inspect centrifuge rotor, buckets, and chamber for damage.																															
Rinse the buckets with lukewarm water and a mild soap or detergent (prior to lubrication).																															
Lubricate the rotor pins and bucket seats with the proper lubricant (e.g., Eppendorf grease).																															
Wipe out the centrifuge chamber with a paper towel sprayed with cleaner.																															
Calibrate the weigh cell.																															
Inspect and clean the waste cabinet.																															
Inspect and clean the reagent cabinet.																															

