

Replacing the air filters

Replace the air filters when an Insufficient hydraulic pressure error is displayed when you attempt to run a sample.

