





The stability of the AMS system is shown in Figure 5, which displays the CCV recoveries during an 8-hour run of 100 simulated seawater samples in Standard mode, with AMS set for 5x dilution. This stability, combined with accurate calibrations at any level, demonstrates the utility of AMS for analysis of samples with challenging matrices.

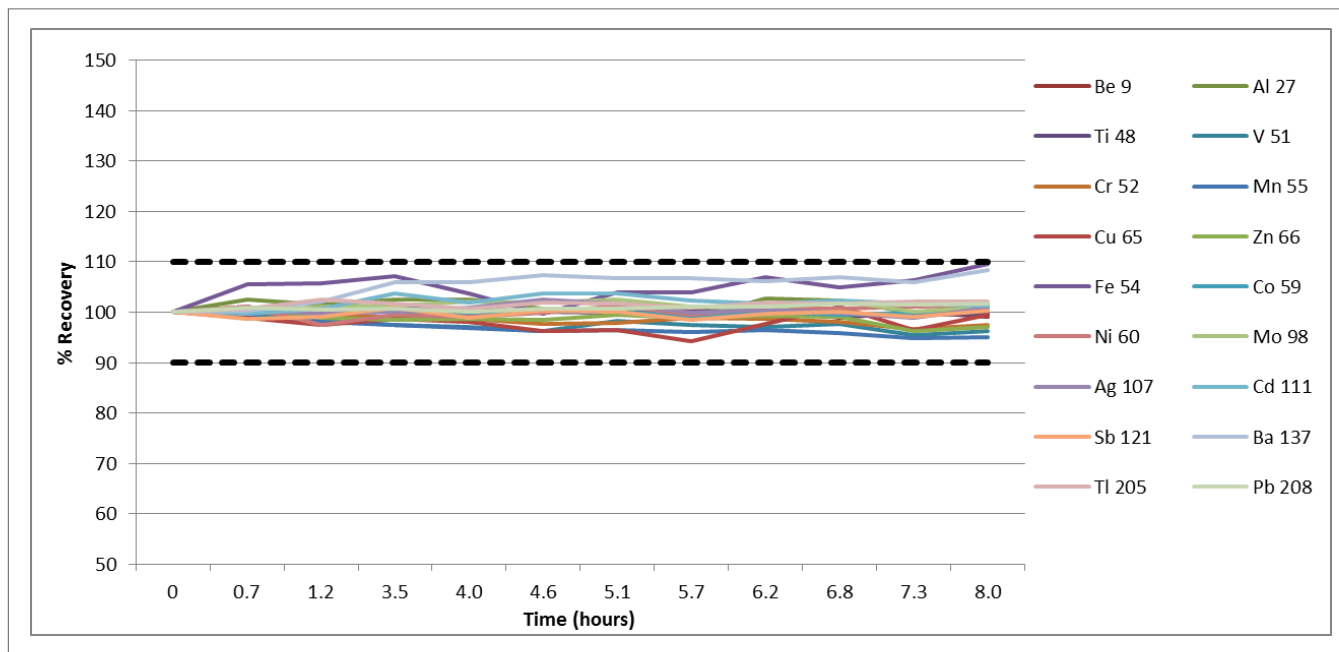


Figure 5. CCV recovery during an 8-hour analysis of 100 simulated seawater samples, with AMS dilution set to 5x.

In summary, PerkinElmer's AMS system provides a number of benefits to simplify analysis of high-matrix samples with the NexION family of ICP-MS instruments. By introducing a flow of argon into the spray chamber neck, the aerosol stream is diluted, allowing for more efficient ionization, fewer matrix effects, and less deposition on the interface cones, which results in simplified sample preparation and higher quality data.