

QCI-706A ULTRAcHECK™ Metals #1 Sample

Please read these instructions carefully before using the quality check sample concentrate

1. Storing the Sample

See package label for information on storage conditions and expiration.

2. Preparing the Sample for Use

Open the vial and using a pipet, transfer 10 mL of the concentrate into a 1 liter volumetric flask (class A) containing approximately 750 mL of reagent water. Add 1 mL concentrated nitric acid, then dilute to volume with reagent water. Mix well. Trace metal grade nitric acid is required. After diluting, use normal holding time and storage conditions for similar samples.

3. Analyzing the Sample

A blind check sample is used to evaluate the quality of the analytical data generated by the laboratory, so use the analytical method your laboratory routinely uses to analyze for these particular analytes.

Samples, standards, and reagent water blanks should all have the same acid matrix. The sample prepared according to the above instructions will contain 0.15% v/v nitric acid; additional or different acids may be used as appropriate for analytical procedures or to match standards.

4. Applicable Methods

Analyte	US EPA Method	Analyte	US EPA Method	Analyte	US EPA Method
Antimony	200.7	Cadmium	200.7	Nickel	200.7
Arsenic	200.7	Chromium	200.7	Selenium	200.7
Barium	200.7	Copper	200.7	Thallium	200.7
Beryllium	200.7	Lead	200.7		

5. Analyte Concentrations

The certificate showing the reference values and advisory ranges is sealed in an envelope, to be opened after the analysis is completed. The advisory ranges represent QC acceptance criteria for analyte recovery following applicable US EPA methodologies. These ranges are based on interlaboratory data, and are included solely as guides for acceptable performance. Each laboratory should develop criteria for judging acceptable method performance based on the intended use of data.