

QCI-702 ULTRAcHECK™ Cations Quality Check 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

Break the ampule, and transfer 10 mL into a 1 liter volumetric flask. Add 20 mL of concentrated nitric acid, and dilute to mark with 18 megohm water. After this product is opened, it should be used and properly disposed of within 2 months. If flame AA is to be used, a further dilution of the above preparation may be necessary. Transfer 5 mL of the above preparation into a 200 mL volumetric flask. Add 2 mL concentrated HNO₃ and dilute to volume with 18 megohm water.

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 method you would normally use to analyze for these particular analytes.

4. Applicable Methods

Analyte	US EPA Method(s)
Calcium	200.7, 215.1
Magnesium	200.7, 242.1
Potassium	200.7, 258.1
Sodium	200.7, 273.1

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.