

Instructions for Use

QCI-750 ULTRAcHECK™ Cyanide 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 open the ampule and transfer 10 mL into a 2 liter glass volumetric flask. Dilute to mark with 18 megohm water. After diluting, this product should be used within 24 hours.

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)
Total Cyanide	335.2, 335.3

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.