

Instructions for Use

QCM-205 ULTRAccheck™ WP Pesticides 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

Add 1000 mL of reagent grade water to a 2 liter separatory funnel. Open the ampule, and transfer 1 mL of the concentrate into the separatory funnel. Mix well. Analyze the sample immediately.

3. Analyzing the Sample

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

4. Applicable Methods

US EPA Method(s) 608, 625, 8080

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