



QCM-251 ULTRAcHECK™ WS Carbamate Pesticides Quality Check Sample

1. Storing the Sample

2. Preparing the Sample for Use

Add approximately 90 mL of reagent water into a 100 mL volumetric flask (class A). Open the ampoule and transfer 1 mL of the concentrate into the flask. Dilute to volume with reagent water. Mix well. Analyze the sample as soon as possible.

3. Analyzing the Sample

4. Applicable Methods

5. Analyte Concentrations

The certificate showing the true values and advisory ranges is sealed in an envelope, to be opened after the analysis is completed. The true values shown for each analyte are the concentrations determined from the gravimetric and volumetric measurements performed during the manufacture of the sample. 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 performance based on the intended use of data.