

# Instructions for Use

## QCM-211      ULTRAccheck™ WS Pesticides #1 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**

The sample concentrate in the ampule must be diluted to its working concentration prior to analysis. Enough sample concentrate is provided to prepare two batches of working sample. Work quickly once the ampule has been opened, as the sample will begin to degrade (through evaporation, for example) once the seal has been broken.

Equilibrate the sample concentrate to 20°C before opening the ampule. Using a syringe or volumetric pipet to transfer the material, add 1.0 mL of the concentrate to 1000 mL of reagent water. Be sure to keep the tip of the syringe 1 cm below the surface of the water while making the addition. Thoroughly mix the working sample prior to use.

### 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.

### 4.      **Applicable Methods**

US EPA Method(s)                      505, 507, 508.1, 525.2

### 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.