

Instructions for Use

QCM-243 ULTRAcHECK™ Aroclors in Transformer Oil Check Sample

Please read these instructions carefully before using the quality check sample.

1. Storing the Sample

See package label for information on storage conditions and expiration.

2. Preparing the Sample for Use

This ULTRAcHECK sample is ready to use. Carefully open the ampule and transfer the required amount of sample.

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. The value listed on the certificate represents the gravimetric true value.

4. Analyte Concentrations

The certificate showing the true values and advisory range 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. The range is based on interlaboratory data, and is included solely as a guide for acceptable performance. Each laboratory should develop criteria for judging acceptable performance based on the intended use of data.