



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY
CINCINNATI, OHIO 45268

Richard F. Jack, Ph.D.
Corporate Market Development
Dionex Corporation
500 Mercury Way
Sunnyvale, CA 94088
408-481-4227

8/15/2008

RE: ATP Case No. D07-0012

Per the terms of the Alternate Test Procedure (ATP) program, the Office of Ground Water and Drinking Water's Technical Support Center (OGWDW/TSC) has determined that the use of chromatographic systems that employ eluent regeneration are acceptable for use in EPA methods 300.0 and 300.1 when operated according to the manufacturer's instructions. Versions of these methods using eluent regeneration systems are determined to be acceptable versions of the original methods as long as all original method quality control parameters are met. Subsequently, these methods may be used for drinking water compliance monitoring performed under National Primary Drinking Water Regulations (NPDWR).

We appreciate your interest in the development of environmental monitoring methods. If you have any questions regarding the review of this alternate test procedure (ATP Case No.D07-0012), please contact Steve Wendelken by e-mail at: wendelken.steve@epa.gov or by telephone at: 513-569-7491.

Sincerely,

A handwritten signature in cursive script that reads "Steve Wendelken".

Steven C. Wendelken PhD.
U.S. EPA, OGWDW-TSC
26 W. Martin Luther King Dr.
Cincinnati, Ohio 45219
Phone: (513) 569-7491
Fax: (513) 569-7837
wendelken.steve@epa.gov

cc:
ATP Coordinators (all Regions)
Danielle Carter, CSC, SCC