



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY  
CINCINNATI, OHIO 45268

Richard F. Jack, Ph.D.  
Corporate Market Development  
Environmental  
Dionex Corporation  
500 Mercury Way  
Sunnyvale, CA  
94088  
[www.dionex.com](http://www.dionex.com)  
408-481-4227

RE: ATP Case Nos. D08-0001, D08-0002

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 electrolytically generated hydroxide eluents and associated hydroxide-selective analytical columns are acceptable for use in EPA methods 317.0 and 326.0. Versions of these method using electrolytically generated hydroxide eluents and associated hydroxide-selective analytical columns 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.D08-0001, D08-002), please contact Steve Wendelken by e-mail at: [wendelken.steve@epa.gov](mailto:wendelken.steve@epa.gov) or by telephone at: 513-569-7491.

Sincerely,

A handwritten signature in cursive script, appearing to read "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](mailto:wendelken.steve@epa.gov)

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