



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY  
CINCINNATI, OHIO 45268

Andrew Eaton, Ph.D.  
Vice President/Laboratory Director  
MWH Laboratories  
750 Royal Oaks Drive, Suite 100  
Monrovia, CA 91016-3629

11/21/2007

RE: ATP Case No. D07-0011

Dear Dr. Eaton

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 MWH analytical methods ML300.0, ML300.1 and ML317.0 [revision dates 11/20/07] are acceptable versions of EPA Methods 300.0, 300.1, and 317.0 respectively. Accordingly, MWH analytical methods ML300.0, ML300.1 and ML317.0 [revision dates 11/20/07] may be used for compliance monitoring performed under National Primary Drinking Water Regulations (NPDWR). MWH analytical methods ML300.0, ML300.1 and ML317.0 [revision dates 11/20/07] add the option of using a quadratic calibration curve to the reference method.

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.D07-0011), 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 black ink, appearing to read "Steve Wendelken".

Steven C. Wendelken, Ph.D.  
ATP Coordinator  
Technical Support Center (MS-140)  
Office of Ground Water and  
Drinking Water

cc:  
ATP Coordinators (all Regions)  
Quality Assurance Managers (all Regions)  
Gregory J. Carroll, USEPA, OGWDW  
Danielle Carter, CSC, SCC