



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY  
WASHINGTON, D.C. 20460

Martin S. Frant, Ph.D.  
Director, Technical Marketing  
Orion Research, Inc.  
500 Cummings Center  
Beverly, MA 01915

MAR - 4 2003

OFFICE OF  
WATER

Dear Dr. Frant:

We are pleased to inform you that the Statistics and Analytical Support Branch (SASB), in concert with the Office of Ground Water and Drinking Water's Technical Support Center (OGWDW/TSC), has determined that Thermo Orion Method AC2009 [Revision 5, April 24, 2002] for determination of fluoride in drinking water (ATP case No. D00-0014) is an acceptable version of EPA-approved methods listed at Title 40 of the *Code of Federal Regulations* (CFR) Part 141 for determining fluoride in drinking water under national primary drinking water regulations (NPDWR).

Method AC2009 determines fluoride after sample distillation using a SPADNS-based colorimetry procedure equivalent to those used in EPA-approved methods, such as Standard Method 4500-F<sup>-</sup>D [20<sup>th</sup> Edition]. Accordingly, Method AC2009 may be used in place of other NPDWR-approved SPADNS-based colorimetry methods for determination of fluoride.

We appreciate your interest in the development of environmental monitoring methods. If you have any questions regarding review of this ATP, please contact Khouane Ditthavong of SASB (202/566-1068) or Herb Brass of OGWDW/TSC (513/569-7936)

Sincerely,

William A. Telliard  
Director of Analytical Methods  
Engineering and Analysis Division (4303)

Herb Brass, Ph.D.  
Technical Support Center (MS-140)  
Office of Ground Water and Drinking Water

cc:

USEPA Regional Administrators (all Regions)  
Quality Assurance Managers (all Regions)  
ATP Coordinators (all Regions)  
Water Management Division Directors (all Regions)  
Gregory J. Carroll, USEPA, OGWDW  
Lillian Holmes, USEPA, OGWDW/TSC  
Maria Gomez-Taylor, USEPA, EAD  
Khouane Ditthavong, USEPA, EAD  
James Boiani, DynCorp, SCC