



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

JUN - 7 2004

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

OFFICE OF  
WATER

Dear Dr. Frant:

We are pleased to inform you that the Statistics and Analytical Support Branch (SASB) has determined that Thermo Orion Method AC2009 [Revision 5, April 24, 2002] for determination of fluoride (ATP case No. N00-0038) is an acceptable version of EPA-approved methods listed at Title 40 of the *Code of Federal Regulations* (CFR) Part 136.3 for determining fluoride in wastewater for National Pollutant Discharge Elimination System (NPDES) compliance monitoring under the Clean Water Act.

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 [18<sup>th</sup>, 19<sup>th</sup>, and 20<sup>th</sup> Editions]. Accordingly, Method AC2009 may be used in place of other NPDES-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 me at 202/566-1061 at your convenience.

Sincerely,

A handwritten signature in black ink, appearing to read "W. A. Telliard".

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

cc:  
USEPA Regional Administrators (all Regions)  
Quality Assurance Managers (all Regions)  
ATP Coordinators (all Regions)  
Water Management Division Directors (all Regions)  
Robin K. Oshiro, USEPA, EAD  
James Boiani, CSC, SCC