



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

JUN - 7 2004

OFFICE OF
WATER

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

Dear Dr. Frant:

We are pleased to inform you that the Statistics and Analytical Support Branch (SASB), has determined that Thermo Orion Method AC2095 [Revision 6, June 4, 2003] for determination of orthophosphate (ATP case No. N00-0040) is an acceptable version of EPA-approved methods listed at Title 40 of the *Code of Federal Regulations* (CFR) Part 136 for determining orthophosphate in wastewater for National Pollutant Discharge Elimination System (NPDES) compliance monitoring under the Clean Water Act.

Method AC2095 determines orthophosphate by colorimetry. Orthophosphate in a sample reacts with ammonium molybdate and potassium antimonyl tartrate in an acid medium to form an antimony-phospho-molybdate complex. The complex is then reduced with ascorbic acid to form intensely colored molybdenum blue which is measured using a colorimeter. The procedure of Method AC2095 is equivalent to those used in EPA-approved methods, such as Standard Method 4500-P E [18th, 19th, and 20th Edition]. Accordingly, Method AC2095 may be used in place of other NPDES-approved colorimetric methods for determination of orthophosphate.

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