



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 AC2046 [Revision 5, April 24, 2002] for determination of nitrite (ATP case No. N00-0039) is an acceptable version of EPA-approved methods listed at Title 40 of the *Code of Federal Regulations* (CFR) Part 136.3 for determining nitrite in wastewater for National Pollutant Discharge Elimination System (NPDES) compliance monitoring under the Clean Water Act.

Method AC2046 determines nitrite by reaction with sulfanilamide to form a diazonium salt which is coupled with N-(1-naphthyl)-ethylenediamine dihydrochloride to form a highly colored azo dye followed by colorimetric detection using a procedure equivalent to those used in EPA-approved methods, such as Standard Method 4500-NO₂-B [18th, 19th, and 20th Edition]. Accordingly, Method AC2046 may be used in place of other NPDES-approved colorimetric methods for determination of nitrite.

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