



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

JUL 24 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 AC2046 [Revision 5, April 24, 2002] for determination of nitrite in drinking water (ATP case No. D00-0015) is an acceptable version of EPA-approved methods listed at Title 40 of the *Code of Federal Regulations* (CFR) Part 141 for determining nitrite in drinking water under national primary drinking water regulations (NPDWR).

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 NPDWR-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 William Telliard of SASB (202/566-1061) or Herb Brass of OGWDW/TSC (513/569-7936) at your convenience.

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
Robin K. Oshiro, USEPA, EAD
James Boiani, DynCorp, SCC