

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 AC4P71 [Revision 4, July 9, 2002] for determination of free chlorine (ATP case No. D02-0004) and Thermo Orion Method AC4P72 [Revision 3, July 9, 2002] for determination of total chlorine (ATP case No. D02-0003) are acceptable versions of EPA-approved methods listed at Title 40 of the Code of Federal Regulations (CFR) Part 141 for national primary drinking water regulations (NPDWR) compliance monitoring.

Methods AC4P71 and AC4P72 determine free chlorine and total chlorine, respectively, using a DPD-based colorimetry procedures equivalent to those in EPA-approved methods such as Standard Method 4500-Cl G [20th Edition]. Therefore Methods AC4P71 and AC4P72 may be used in place of NPDWR-approved DPD-based colorimetry methods for free and total chlorine, respectively.

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