



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

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

SEP 3 1 2002

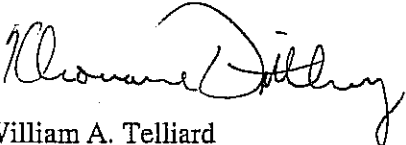
OFFICE OF
WATER

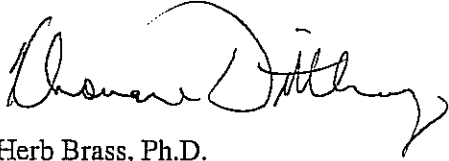
Dear Dr. Frant:

We are pleased to inform you that the Statistics and Analytical Support Branch (SASB) and the Office of Ground Water and Drinking Water's Technical Support Center (OGWDW/TSC) has determined that Thermo Orion Method AC2072: "Low Range Total Chlorine" [March 14, 2002, Revision 4] (ATP Case No. D00-0013) is an acceptable version of EPA-approved methods listed at Title 40 of the *Code of Federal Regulations* (CFR) Part 141 for determining total residual chlorine in drinking water by reaction with N, N-diethyl-p-phenyldiamine sulfate (DPD) followed by spectrophotometric detection (e.g., Standard Method 4500-Cl G). Accordingly, Thermo Orion Method AC2072 may be used for NPDWR compliance monitoring. Both the approved methods and the Thermo Orion method rely on the same chemistry (reaction of oxidants with DPD) to produce a red color; the intensity of which is: (1) dependent upon chlorine concentration and (2) determined photometrically.

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,


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


for 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