

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

MAR 20 2002

OFFICE OF

Dr. Martin S. Frant 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 AC2072: "Low Range Total Chlorine" [March 14, 2002, Revision 4] (ATP Case No. N00-0037) is an acceptable version of EPA-approved methods listed at Title 40 of the Code of Federal Regulations (CFR) Part 136.3, Table IB, for determining total residual chlorine 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 NPDES compliance monitoring. Both the approved methods and the Thermo Orion method rely on the same chemistry (i.e., 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 (202/260-6115).

Sincerely,

William A. Telliard, Director

Analytical Methods

Engineering and Analysis Division (4303)

ce:

Maria Gomez-Taylor, USEPA, EAD
Khouane Ditthavong, USEPA, EAD
USEPA Regional Administrators (all Regions)
Quality Assurance Managers (all Regions)
ATP Coordinators (all Regions)
Water Management Division Directors (all Regions)