



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY

WASHINGTON, D.C. 20460

NOV 19 2002

Douglas W. Later, Ph.D.
Environmental Market Development Manager
Dionex Corporation
500 Mercury Drive
Sunnyvale, California 94086

OFFICE OF
WATER

Dear Dr. Later:

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) have determined that the use of hydroxide eluents (either manually prepared or electrolytically generated) falls within the method flexibility allowed in EPA Methods 300.0 and 300.1 for determining inorganic anions. Specifically, Sections 9.4.6 and 11.8 of Method 300.0 (Rev. 2.1, August 1993) and Section 9.4.5 of Method 300.1 (Rev. 1.0, 1997) state that the substitution of eluents to improve method performance is an acceptable method modification. Therefore, the use of hydroxide eluents in EPA Methods 300.0 and 300.1 is acceptable for compliance monitoring under the Clean Water Act and Safe Drinking Water Act. This determination supersedes all prior determinations on this subject and applies to all revisions of EPA Methods 300.0 and 300.1.

If you have any questions regarding this determination, 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 (4303T)

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
James Boiani, DynCorp, SCC
Robert J. Joyce, Dionex