



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY

WASHINGTON, D.C. 20460

Wilbur H. (Bill) Campbell
President
The Nitrate Elimination Co., Inc (NECi).
334 Hecla St., Lake Linden MI 49945
Bill@nitrate.com

OFFICE OF WATER

SUBJECT: Review of NECi Method N07-0003, Method for Nitrate Reductase Nitrate-Nitrogen Analysis (ATP Case No. N07-0003)

DATE: April 4, 2014

I have reviewed NECi Method N07-0003 (ATP Case No. N07-003), "Method for Nitrate Reductase Nitrate-Nitrogen Analysis", and the supporting validation data in ATP Case No. N07-0003. I determined that this method meets all requirements for measurement of nitrate and combined nitrate-nitrite in wastewater. That is, the performance of this method is substantially similar to methods listed at 40 CFR Part 136 for measurement of nitrate and combined nitrate-nitrite in wastewater.

I will recommend that this method be included in future regulatory actions in which EPA periodically adds to the list of approved methods at 40 CFR Part 136. However, this ATP review does not replace this notice and comment rulemaking process. In the interim a user may, on a facility-by-facility basis, seek approval from their regional authority for use of this method in measuring nitrate and combined nitrate-nitrite in Clean Water Act (CWA) programs.

If I can be of any additional assistance on this matter or others, please contact me at walker.lemuel@epa.gov.

Sincerely,

A handwritten signature in blue ink that reads "Lemuel Walker, Jr.".

Lemuel Walker, Jr.
CWA ATP Coordinator
Engineering and Analytical Support Branch
Engineering and Analysis Division
Office of Science and Technology

cc: Quality Assurance Managers (all Regions)
ATP Coordinators (all Regions)