



UNITED STATES ENVIRONMENTAL PROTECTION AGENCY
CINCINNATI, OHIO 45268

Richard F. Jack, Ph.D.
Corporate Market Development
Environmental
Dionex Corporation
500 Mercury Way
Sunnyvale, CA
94088
www.dionex.com
408-481-4227

Richard

To provide clarification, capillary chromatography columns are an acceptable alternative for any EPA method that allows flexibility in column choice. Although Method 300.1 gives the example of 2mm or 4mm in section 6.1.2 capillary columns are acceptable as long as all of the quality control parameters of the method are met. This flexibility is already built into the methods and does not require additional approval.

We appreciate your interest in the development of environmental monitoring methods. If you have any questions regarding this issue please contact Steve Wendelken by e-mail at: wendelken.steve@epa.gov or by telephone at: 513-569-7491.

Sincerely,

A handwritten signature in black ink, appearing to read "Steve Wendelken".

Steven C. Wendelken PhD.
U.S. EPA, OGWDW-TSC
26 W. Martin Luther King Dr.
Cincinnati, Ohio 45219
Phone: (513) 569-7491
Fax: (513) 569-7837
wendelken.steve@epa.gov

cc:
ATP Coordinators (all Regions)
Danielle Carter, CSC, SCC