E. Kevin Hrusovsky Chairman, President and Chief Executive Officer Quanterix Corporation 113 Hartwell Avenue Lexington, MA 02421

> Re: Quanterix Corporation Registration Statement on Form S-3 File No. 333-230399 Filed March 19, 2019

Dear Mr. Hrusovsky:

This is to advise you that we have not reviewed and will not review your registration statement.

Please refer to Rules 460 and 461 regarding requests for acceleration. We remind you that the company and its management are responsible for the accuracy and adequacy of their disclosures, notwithstanding any review, comments, action or absence of action by the staff.

Please contact Thomas Jones at 202-551-3602 with any questions.

Sincerely, Division of Office of

Corporation Finance

Electronics and Machinery cc: William T. Whelan, Esq.