

UNITED STATES
SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

Notice of Effectiveness

Effectiveness Date: May 10, 2019 5:30 P.M.

Form: S-3

CIK: [0001503274](#)

Company Name: Quanterix Corp

File Number: [333-230399](#)