

UNITED STATES
SECURITIES AND EXCHANGE COMMISSION
Washington, D.C. 20549
Notice of Effectiveness

Effectiveness Date: June 12, 2025
Accession Number: 0001104659-25-056131
Submission Type: POS AM

CIK: [0001503274](#)
Company Name: Quanterix Corp
File Number: [333-284932](#)