## UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

Notice of Effectiveness	
<b>Effectiveness Date:</b>	December 6, 2017 5:05 P.M.
Form:	S-1
CIK:	0001503274
Company Name:	Quanterix Corp
File Number:	<u>333-221475</u>
Ĩ	