


Product Number:	101736 (103159)
Lot Number:	501554 (83410)
Expiration:	7-Dec-2019
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval Sivasankari Murugan QC Supervisor  14DEC2018

Name Title Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.