

Product Number:	101736 (103159)
Lot Number:	501417 (82350)
Expiration:	23-Aug-2019
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass


Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title


Signature/Date

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