

Product Number:	101736 (103159)
Lot Number:	501237 (81022)
Expiration:	12-Apr-2019
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 02 MAY 2018
Signature/Date

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