


Product Number:	101736 (103159)
Lot Number:	501477 (82910)
Expiration:	18-Oct-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.