

Product Number:	101736 (103159)
Lot Number:	501317 (81730)
Expiration:	22-Jun-2019
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

  
Signature/Date

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