

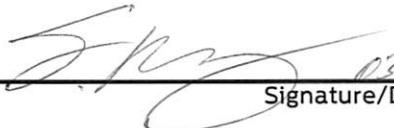
|                    |                 |
|--------------------|-----------------|
| Product Number:    | 101736 (103159) |
| Lot Number:        | 501357 (81930)  |
| Expiration:        | 12-Jul-2019     |
| Storage Condition: | 2-8°C           |

|                  | PASS/FAIL |
|------------------|-----------|
| Release Criteria | Pass      |

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

 03 Aug 2018  
Signature/Date

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