

<b>Product Number:</b>	101736 (103159)
<b>Lot Number:</b>	501397 (81650)
<b>Expiration:</b>	14-Jun-2019
<b>Storage Condition:</b>	2-8°C

	<b>PASS/FAIL</b>
<b>Release Criteria</b>	Pass

**Review/Approval**

Sivasankari Murugan  
**Name**

QC Supervisor  
**Title**

  
**Signature/Date** 09 SEP 2018

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