

<b>Product Number:</b>	101736 (103159)
<b>Lot Number:</b>	501701 (90600)
<b>Expiration:</b>	1-Mar-2020
<b>Storage Condition:</b>	2-8°C

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

Review/Approval

Patrizia Stadler

Associate Director, QC

*p. stadler 07 Mar 2019*

Name

Title

Signature/Date

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