

<b>Product Number:</b>	101736 (100030)
<b>Lot Number:</b>	500731 (70701)
<b>Expiration:</b>	10-Dec-2017
<b>Storage:</b>	2-8°C

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

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler*

19 Sep 2019

Name

Title

Signature/Date