

<b>Product Number:</b>	101736 (103159)
<b>Lot Number:</b>	502232 (00170)
<b>Expiration:</b>	17-Jan-2021
<b>Storage:</b>	2-8°C

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

Review/Approval

Patrizia Stadler Assoc. Director, QC  
Name Title

*P Stadler*

10 Feb 2020

Signature/Date