

<b>Product Number:</b>	101736 (103159)
<b>Lot Number:</b>	502199 (93460)
<b>Expiration:</b>	12-Dec-2020
<b>Storage:</b>	2-8°C

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

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	14 Jan 2020
<b>Name</b>	<b>Title</b>	<b>Signature</b>	<b>Date</b>