

Product Number:	101736 (103159)
Lot Number:	502151 (93090)
Expiration:	5-Nov-2020
Storage:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	16 Dec 2019
Name	Title	Signature	Date