

Product Number:	101736 (103159)
Lot Number:	502260 (00290)
Expiration:	29-Jan-2021
Storage:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

28 Feb 2020

Name

Title

Signature/Date