

Product Number:	100033
Lot Number:	104601
Expiration:	15-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval

Ruby Burke
Name

QC, Sr. Director
Title

Ruby Burke 05mar2021
Signature/Date