

Product Number:	103397
Lot Number:	104620
Expiration:	15-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 11 Mar 2021
Signature/Date