

Product Number:	103397
Lot Number:	120031
Expiration:	28-Jun-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivam Goel
Name

QC
Title

Shivam Goel 17 Aug 2021
Signature/Date