

Product Number:	103397
Lot Number:	116813
Expiration:	17-Jun-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 14 Jul 2021
Signature/Date