

Product Number:	103650
Lot Number:	106718
Expiration:	8-Jun-2021
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 22Mar2021
Signature/Date