

Product Number:	103650
Lot Number:	127010
Expiration:	27-Dec-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	139
---------------------------------------	-----

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 13Oct2021
Signature/Date