

Product Number:	103397
Lot Number:	301009
Expiration:	10-Jan-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Sonia Khilka
Name

QC Manager
Title

Sonia Khilka 06 APR 2023
Signature/Date