

Product Number:	101735
Lot Number:	503540
Expiration:	1-Jun-2023



Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	215209
SBG Concentrate	103397	215710

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Della / 08 JUL 2022
Signature/Date