

Product Number:	101735
Lot Number:	503804
Expiration:	29-Aug-2023



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

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	302512
SBG Concentrate	103397	230622

Laboratory Analysis

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

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 08 FEB 2023
Signature/Date