

Product Number:	101735
Lot Number:	503392
Expiration:	19-Oct-2022



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

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	203113
SBG Concentrate	103397	129210

Laboratory Analysis

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 08 Mar 2022
Signature/Date