

Product Number:	104074
Lot Number:	503700
Paired Reagent Kit	503501
Expiration:	1-May-2023
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	104071	218805
Calibrator Diluent	104072	212103
Calibrator Concentrate	102619	220014

Calibrator Concentrate Value Assignment (pg/mL)	3,887
---	-------



Data below represents results generated on the Simoa[®] HD-X Analyzer.

Review/Approval

Sonia Khillan

QC

Sonia Khillan / 130CT2022