

Product Number:	104074
Lot Number:	503974
Paired Reagent Kit	503808
Expiration:	31-Jan-2024
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	104071	304005
Calibrator Diluent	104072	303110
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 Khilka

QC

Sonia Khilka 22SEP2023