

Product Number:	103715
Lot Number:	502616
Paired Reagent Kit	502613
Expiration:	13-May-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	103704	031808
Calibrator Diluent	103707	031520
Calibrator Concentrate	103705	027201

Calibrator Concentrate Value Assignment (pg/mL)	1223
--	------

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 03 Dec 20 20
Signature/Date