

Product Number:	103807
Lot Number:	502709
Paired Reagent Kit	502708 (Product No. 103806)
Expiration:	25-Jan-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	103800	102524
Calibrator Diluent	103802	102526
Calibrator Concentrate	103801	105602

Calibrator Concentrate Value Assignment (pg/mL)	3840
---	------

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 01 Mar 2021
Signature/Date