

Product Number:	103715
Lot Number:	502558
Paired Reagent Kit	502539
Expiration:	18-Mar-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Sample Diluent	103704	026201
Calibrator Diluent	103707	026722
Calibrator Concentrate	103705	027201

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

Review/Approval

Shivani Goel

QC Manager

Name

Title

Shivani Goel 06Oct2020

Signature/Date