

Product Number:	103346
Lot Number:	5330120250
Expiration Date:	23-Aug-2027

Component	Item Number	Lot Number
Simoa® Cuvettes	103160	E24083G6

Characteristic Inspected	PASS / FAIL
Raw Material	PASS
Visual	PASS
Physical	PASS

Review/Approval

Dipesh Podd
Name

Sr. QA Specialist
Title

DT 21Aug2025
Signature/Date