

**CERTIFICATE OF ANALYSIS**

Product Description: S 600/600MM and SQ 120/120MM PM/Installation Kit
Catalog Number: RA1IM-1
Lot Number: PI100098
Expiration Date: 30 JUN 2027
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
96-Well 1-Spot SECTOR Plate	21Q1EA	31 MAR 2028
MULTI-SPOT® 96-well 10-spot X-talk Plate	2BPY2A	31 OCT 2029
96-well 10-Spot Plate Assy, STD, SECTOR	2BPGPA	30 JUN 2027
384-well 1-Spot Plate Assy, STD, SECTOR	05P0DA	31 JUN 2027
MSD Free Tag ECL 300,000	Y0250051	30 SEP 2028
MSD Free Tag ECL 15,000	Y0260187	28 FEB 2029
PQ Control 2	Y0810011	30 NOV 2028

MSD Testing Results:

Metric	Observed
MSD Free Tag 15000 Mean Signal	15035

Specifications:

Specification	Minimum	Maximum
MSD Free Tag 15000 Mean Signal	12000	18000
MSD Free Tag 15000 %CV	0	6.0%

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		27 JUL 2026