



CERTIFICATE OF ANALYSIS

Product Description: S 600/600MM and SQ 120/120MM PM/Installation Kit
Catalog Number: RA1IM-1
Lot Number: PI100097
Expiration Date: 30 JUN 2027
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
96-Well 1-Spot SECTOR Plates	21Q0EA	31 JAN 2028
MULTI-SPOT® 96-well 10-spot X-talk Plate	2BPY2A	31 OCT 2029
MULTI-SPOT® 96-well 10-spot HB Plate	2CQ01A	31 JAN 2028
MULTI-ARRAY® 384-well 1-spot Plate	05P0TA	30 JUN 2027
MSD Free Tag ECL 300,000	Y0250051	30 SEP 2028
MSD Free Tag ECL 15,000	Y0260187	28 FEB 2029

MSD Testing Results:

Metric	Observed
MSD Free Tag 15000 Mean Signal	14538

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	Karen Hamilla	Quality		05 JUN 2026