



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S 600/600MM and SQ 120/120MM PM/Installation Kit

Catalog Number: RA1IM-1

Lot Number: PI100078

Expiration Date: 28 FEB 2025

Storage Temperature: Room Temperature

Specifications:

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

MSD Testing Results:

Metric	Observed
MSD Free Tag 15000 Mean Signal	14058

Lot-specific Components:

Description	Lot Number	Expiration Date
MULTI-ARRAY® 96-well 1-spot plate	21N20A	30 APR 2025
MULTI-SPOT® 96-well 10 spot X-talk Plate	2BNY1A	28 FEB 2027
MULTI-SPOT 96-well 10 spot HB Plate	2CN0QA	30 JUN 2025
MULTI-ARRAY 384-well 1-spot Plate	05N02A	28 FEB 2025
MSD Free Tag ECL 15000	Y0260178	28 FEB 2026
MSD Free Tag ECL 300000	Y0250046	31 DEC 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 AUG 2023