



CERTIFICATE OF ANALYSIS

Product Description: S 600/600MM and SQ 120/120MM PM/Installation Kit
Catalog Number: RA1IM-1
Lot Number: PI100080
Expiration Date: 28 FEB 2025
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
MULTI-ARRAY® 384-well 1-spot Plate	05N02A	28 FEB 2025
MULTI-ARRAY® 96-well 1-spot plate	21N2EA	31 MAY 2025
MULTI-SPOT® 96-well 10-spot HB Plate	2CN0SA	30 JUN 2025
MULTI-SPOT® 96-well 10-spot X-talk Plate	2BNY2A	30 APR 2027
MSD Free Tag ECL 300,000	Y0250047	31 MAY 2026
MSD Free Tag ECL 15,000, 250mL	Y0260179	31 AUG 2026

MSD Testing Results:

Metric	Observed
MSD Free Tag 15000 Mean Signal	14280

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		15 DEC 2023