



### 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:** PI100079

**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	14280

**Lot-specific Components:**

Description	Lot Number	Expiration Date
MULTI-ARRAY® 96-well 1-spot plate	21N2EA	31 MAY 2025
MULTI-SPOT® 96-well 10 spot X-talk Plate	2BNY2A	30 APR 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	Y0260179	31 AUG 2026
MSD Free Tag ECL 300000	Y0250047	31 MAY 2026

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		18 OCT 2023