

**CERTIFICATE OF ANALYSIS**

**Product Description:** S 600/600MM and SQ 120/120MM PM/Installation Kit  
**Catalog Number:** RA1IM-1  
**Lot Number:** PI100082  
**Expiration Date:** 30 JUN 2025  
**Storage Temperature:** Room Temperature

**Lot-Specific Components:**

Description	Lot Number	Expiration Date
MULTI-ARRAY® 384-well 1-spot Plate	05N0JA	30 JUN 2025
MULTI-ARRAY® 96-well 1-spot plate	21N3MA	31 JUL 2025
MULTI-SPOT® 96-well 10-spot HB Plate	2CN0VA	30 JUN 2025
MULTI-SPOT® 96-well 10-spot X-talk Plate	2BNY3A	30 JUN 2027
MSD Free Tag ECL 300,000	Y0250047	31 MAY 2026
MSD Free Tag ECL 15,000, 250mL	Y0260180	31 DEC 2026

**MSD Testing Results:**

Metric	Observed
MSD Free Tag 15000 Mean Signal	14631

**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
<b>Review/Approval</b>	Karen Hamilla	Quality		14 MAR 2024