



CERTIFICATE OF ANALYSIS

**Product Description:** S 600/600MM and SQ 120/120MM PM/Installation Kit  
**Catalog Number:** RA1IM-1  
**Lot Number:** PI100096  
**Expiration Date:** 30 JUN 2027  
**Storage Temperature:** Room Temperature

**Lot-Specific Components:**

| Description                              | Lot Number | Expiration Date |
|--|------------|-----------------|
| 96-Well 1-Spot SECTOR Plates             | 21P50A     | 31 OCT 2027     |
| MULTI-SPOT® 96-well 10-spot X-talk Plate | 2BOY2A     | 30 NOV 2028     |
| MULTI-SPOT® 96-well 10-spot HB Plate     | 2CQ01A     | 31 JAN 2028     |
| MULTI-ARRAY® 384-well 1-spot Plate       | 05P0LA     | 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 | 14565    |

**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 MAY 2026 |