

**CERTIFICATE OF ANALYSIS**

Product Description: S 600/600MM and SQ 120/120MM Performance Qualification Kit
Catalog Number: R31QQ-3
Lot Number: PQ300160
Expiration Date: 31 MAY 2025
Storage Temperature: Room Temperature

Lot-Specific Components:

| Description | Lot Number | Expiration Date |
|----------------------------|------------|-----------------|
| MULTI-ARRAY® 96-Well Plate | 21N2EA | 31 MAY 2025 |
| MSD Free Tag ECL 15000 | Y0260179 | 31 AUG 2026 |
| PQ Low Control | Y0110099 | 31 JUL 2026 |

MSD Testing Results:

| Metric | Observed |
|----------------------------|----------|
| Free Tag Mean Signal | 14280 |
| PQ Low Control Mean Signal | 69 |
| Dark Noise Mean Signal | 3 |

Specifications:

| Specification | Minimum | Maximum |
|-----------------------------------|---------|---------|
| Free Tag Mean Signal | 12000 | 18000 |
| Free Tag %CV | 0 | 6% |
| PQ Low Control Mean Signal | 25 | 100 |
| PQ Low Control Standard Deviation | 0 | 25 |
| Dark Noise Mean Signal | -16 | 16 |
| Dark Noise Standard Deviation | 0 | 16 |

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

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 17 OCT 2023 |