



CERTIFICATE OF ANALYSIS

**Product Description:** S 600/600MM and SQ 120/120MM Performance Qualification Kit  
**Catalog Number:** R31QQ-3  
**Lot Number:** PQ300163  
**Expiration Date:** 28 FEB 2026  
**Storage Temperature:** Room Temperature

**Lot-Specific Components:**

| Description                | Lot Number | Expiration Date |
|----------------------------|------------|-----------------|
| MULTI-ARRAY® 96-Well Plate | 21O0NA     | 28 FEB 2026     |
| 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       | 14052    |
| PQ Low Control Mean Signal | 57       |
| 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        |
|------------------------|---------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |           | 21 MAR 2024 |