



CERTIFICATE OF ANALYSIS

Product Description: S 600/600MM and SQ 120/120MM Performance Qualification Kit
Catalog Number: R31QQ-3
Lot Number: PQ300183
Expiration Date: 31 OCT 2027
Storage Temperature: Room Temperature

Lot-Specific Components:

| Description | Lot Number | Expiration Date |
|------------------------------|------------|-----------------|
| 96-Well 1-Spot SECTOR Plates | 21P5OA | 31 OCT 2027 |
| MSD Free Tag ECL 15000 | Y0260187 | 28 FEB 2029 |
| PQ Low Control | Y0320018 | 30 SEP 2028 |

MSD Testing Results:

| Metric | Observed |
|----------------------------|----------|
| Free Tag Mean Signal | 14565 |
| PQ Low Control Mean Signal | 54 |
| Dark Noise Mean Signal | 1 |

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 | | 14 MAY 2026 |