



CERTIFICATE OF ANALYSIS: TITRATION PLATES

COA – 8633-B

Product Description: MSD GOLD™ 96-Well Small Spot Streptavidin SECTOR® Plate

Catalog #: L45SA-1 (1 plate); L45SA-2 (5 plates); L45SA-5 (30 plates);
L45SA-6 (120 plates); L45SA-7 (510 plates)

Lot Number: Z0021316

Expiration date: 31 DEC 2019

Statement: This product was manufactured and tested according to approved MSD Standard Operating Procedures.

This product has an approximate binding capacity of 0.075 pmoles of biotinylated IgG. For optimal assay results, the amount of biotinylated reagent used in the assay should not exceed this binding capacity.

Titration plates were tested using tagged BTI lot: R0010578

| Concentration of SULFO-TAG™ Conjugated Biotinylated IgG (BTI) | Mean ECL spec and tolerance range | Pass/Fail? |
|--|-----------------------------------|------------|
| 1.5 nM (0.075 pmole) | 45,472 counts ± 15% | Pass |
| 1.0 nM (0.05 pmole) | 34,498 counts ± 15% | Pass |
| 0.5 nM (0.025 pmole) | 18,826 counts ± 15% | Pass |
| 0 nM (0.0 pmole) | Not to exceed 100 counts | Pass |

| Concentration of SULFO-TAG™ Conjugated Biotinylated IgG (BTI) | Intraplate CV or SD spec | Pass/Fail? |
|--|--------------------------|------------|
| 1.5 nM (0.075 pmole) | ≤10% | Pass |
| 1.0 nM (0.05 pmole) | ≤10% | Pass |
| 0.5 nM (0.025 pmole) | ≤10% | Pass |
| 0 nM (0.0 pmole) | ≤15 count SD | Pass |



CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

COA - 8633-B

This product has an approximate binding capacity of 0.075 pmoles of biotinylated IgG. For optimal assay results, the amount of biotinylated reagent used in the assay should not exceed this binding capacity.

| Metric | Specification | Pass/Fail? |
|---|---|------------|
| Mean Intraplate CV | $\leq 6\%$ | Pass |
| Intraplate CV | $\leq 8\%$ for at least 91.5% of plates | Pass |
| Number of plates with intraplate CV $>12\%$ | 0 plates | Pass |
| Interplate CV | $\leq 8\%$ | Pass |
| Plates where signal $> 20\%$ from plate mean occurs in same well on multiple plates | 0 plates | Pass |
| Wells with signal $> 50\%$ from plate mean | 0 plates | Pass |
| Median signal for concentric rings, min to max range | $\leq 10\%$ for all plates | Pass |
| Median signal for columns, min to max range | $\leq 10\%$ for at least 91.5% of plates | Pass |
| Median signal for columns, min to max range | $\leq 15\%$ for all plates | Pass |
| Median signal for rows, min to max range | $\leq 10\%$ for at least 91.5% of plates | Pass |
| Median signal for rows, min to max range | $\leq 15\%$ for all plates | Pass |
| Number of plates sampled (N) | Lot size dependent; see SOP 3-2000-6P0002 | Pass |

The above product is intended for research use only. Not for use in diagnostic procedures.

| | Name | Function | Signature | Date |
|-----------------|-----------------|-------------------|-----------|-------------|
| Review/Approval | Carissa Hubbard | Quality Control | | 05 OCT 2017 |
| Review/Approval | Dawn Stollar | Quality Assurance | | 05 OCT 2017 |