



CERTIFICATE OF ANALYSIS: TITRATION PLATES

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

Catalog #: L45SA - Series

Lot Number: Z0021600

Date of Manufacture: 12 NOV 2019

Expiration Date: 30 APR 2022

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

This product has an approximate binding capacity of 0.075 pmol 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: R0010605

| Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI) | Referenced Signal | Pass/Fail? |
|--|--------------------------|------------|
| 1.5 nM (0.075 pmol) | 45472 counts ± 15% | Pass |
| 1 nM (0.05 pmol) | 34498 counts ± 15% | Pass |
| 0.5 nM (0.025 pmol) | 18826 counts ± 15% | Pass |
| 0 nM (0.0 pmol) | Not to Exceed 100 Counts | Pass |

| Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI) | Intraplate CV or SD | Pass/Fail? |
|--|---------------------|------------|
| 1.5 nM (0.075 pmol) | ≤ 10% | Pass |
| 1 nM (0.05 pmol) | ≤ 10% | Pass |
| 0.5 nM (0.025 pmol) | ≤ 10% | Pass |
| 0 nM (0.0 pmol) | ≤ 15 Count SD | Pass |



CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

Statement: This product has an approximate binding capacity of 0.05 pmol 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 CVs | $CV \leq 8\%$ for 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 Wells | Pass |
| Median signal for concentric rings, min to max range | $\leq 10.0\%$ for 100% of Plates | Pass |
| Median signal for columns, min to max range | $\leq 10.0\%$ for 91.5% of Plates | Pass |
| Median signal for columns, min to max range | $\leq 15.0\%$ for 100% of Plates | Pass |
| Median signal for rows, min to max range | $\leq 10.0\%$ for 91.5% of Plates | Pass |
| Median signal for rows, min to max range | $\leq 15.0\%$ for 100% of Plates | Pass |

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

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 20 NOV 2019 |