



CERTIFICATE OF ANALYSIS: TITRATION PLATES

Product Description: MSD GOLD™ 96-well Streptavidin SECTOR® Plate

Catalog #: L15SA - Series

Lot Number: Z0021549

Date of Manufacture: 12 AUG 2019

Expiration Date: 31 JAN 2022

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

This product has an approximate binding capacity of 0.3 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? |
|--|--------------------------|------------|
| 6 nM (0.3 pmol) | 57096 counts ± 15% | Pass |
| 4 nM (0.2 pmol) | 46132 counts ± 15% | Pass |
| 2 nM (0.1 pmol) | 25185 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? |
|--|---------------------|------------|
| 6 nM (0.3 pmol) | ≤ 10% | Pass |
| 4 nM (0.2 pmol) | ≤ 10% | Pass |
| 2 nM (0.1 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 0.2 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 | Xiomara Talavera | Quality | | 29 AUG 2019 |