



## CERTIFICATE OF ANALYSIS: TITRATION PLATES

COA - 08008

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

**Catalog #:** L15SA-1 (1 plate); L15SA-2 (5 plate); L15SA-5 (30 plate);  
L15SA-6 (120 plate); L15SA-7 (510 plate)

**Lot Number:** Z0021276

**Expiration date:** 31Jul2019

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

This product has an approximate binding capacity of 0.3 pmole 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™ Labeled Biotinylated IgG (BTI) | Mean ECL spec and tolerance range | Pass/Fail? |
|--|-----------------------------------|------------|
| 6 nM (0.3 pmole)   | 66,097 counts ± 15%               | Pass       |
| 4 nM (0.2 pmole)   | 50,830 counts ± 15%               | Pass       |
| 2 nM (0.1 pmole)   | 25,531 counts ± 15%               | Pass       |
| 0 nM (0.0 pmole)   | Not to exceed 100 counts          | Pass       |

| Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI) | Intraplate CV or SD spec | Pass/Fail? |
|--|--------------------------|------------|
| 6 nM (0.3 pmole)   | 10%                      | Pass       |
| 4 nM (0.2 pmole)   | 10%                      | Pass       |
| 2 nM (0.1 pmole)   | 10%                      | Pass       |
| 0 nM (0.0 pmole)   | 15 count SD              | Pass       |



## CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

COA - 08008

This product has an approximate binding capacity of 0.3 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 CVs  | $\leq 8\%$ 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 | Xiomara Talavera | Quality Control   | <i>Talavera</i>     | 17 MAR 2017 |
| Review/Approval | Dawn Stollar     | Quality Assurance | <i>Dawn Stollar</i> | 17 MAR 2017 |