



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD 96-Well 1-Spot Streptavidin SECTOR Plate
Catalog Number: L15SA-Series
Lot Number: Z0022383
Expiration Date: 30 APR 2027

The above plate lot was tested according to MSD's procedures and passed all specifications. Plates were tested using SULFO-TAG labeled biotinylated IgG (BTI).

Titration Results:

Table with 3 columns: Metric, Specification, Result. Rows include 0.300 pmol BTI, 0.200 pmol BTI, 0.100 pmol BTI, Max Background Signal, and Standard Deviation of Background Signal.

Uniformity Results:

Plates sampled from the above plate lot were tested by adding 0.200 pmol BTI to each well.

Table with 3 columns: Metric, Specification, Result. Rows include Average Intraplate CV, Percentage of plates with CV <= 8.0%, Number of plates with Intraplate CV > 12%, and Interplate CV.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

Table with 5 columns: Review/Approval, Name, Function, Signature, Date. Row includes Karen Hamilla, Quality, signature, and 22 NOV 2024.