



CERTIFICATE OF ANALYSIS

Product Description: MSD Read Buffer T (4x), Surfactant Free
CatalogNumber/Size: R92TD-Series
Lot Number: Y0060032
Expiration Date: 31 Aug 2022
Storage Temperature: Room Temperature

Statement: MSD Read Buffer T (4x), Surfactant Free is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD Read Buffer T (4x), Surfactant Free are provided in the product insert available at www.mesoscale.com.

MSD Read Buffer T (4x) Surfactant Free was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010605 and MSD GOLD™ 96-well High Bind Avidin SECTOR Plate lot # Z0010663

| Concentration of BTI | Mean ECL spec and tolerance range | Intraplate CV or SD spec | Pass/Fail? |
|----------------------|-----------------------------------|--------------------------|------------|
| BTI with 4X | 48609 - 72914 counts | 15 % | PASS |
| Dil-100 with 4X | 0 - 158 counts | 50 % | PASS |

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|----------------------|-------------|
| Review/Approval | Karen Hamilla | Quality | <i>Karen Hamilla</i> | 16 DEC 2019 |