



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD Read Buffer T (4X) With Surfactant

Catalog #/Size: R92TC-Series

Lot Number: Y0140363

Expiration Date: 30 NOV 2021

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Signal test, using controlled reagents and plates:

ECL, positive control (tolerance range): 1X: 25 565 – 47 479
2X: 43 205 – 58 454

Baseline ECL (tolerance range): 1X: 19 – 44
2X: 38 – 88

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

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 10 DEC 2018 |