



CERTIFICATE OF ANALYSIS

COA - 07166

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

Catalog #/Size: R92TC-1 / 1000 mL; R92TC-2 / 200 mL

Lot Number: Y0140336

Expiration Date: 30 JUN 2019

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

pH: 7.7 ± 0.1

Signal test, using controlled reagents and plates:

ECL, positive control (mean \pm tolerance): 1X: $36\,522 \pm 30\%$

2X: $50\,830 \pm 15\%$

Baseline ECL: 1X: $31.5 \pm 40\%$

2X: $63 \pm 40\%$

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

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>Xiomara Talavera</i>	27 JUL 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 JUL 2016