



## CERTIFICATE OF ANALYSIS

COA - 06767

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

**Catalog #/Size:** R92TC-1 / 1000 mL; R92TC-2 / 200 mL; R92TC-3 / 50 mL

**Lot Number:** Y0140330

**Expiration Date:** 31 JAN 2019

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

pH:  $7.8 \pm 0.1$

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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		11 APR 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		11 APR 2016