



CERTIFICATE OF ANALYSIS

COA - 07071

Product Description: MSD Read Buffer T (4x) with Surfactant

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

Lot Number: Y0140333

Expiration Date: 31 MAR 2019

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

pH: 7.8 ± 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	Rolanda Spratley	Biologics/Chemistry		01 JUL 2016
Review/Approval	Dawn Stollar	Quality Assurance		01 JUL 2016