



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD Read Buffer B
Catalog #/Size: R92BA-Series
Lot Number: Y0530001
Expiration Date: 31 DEC 2020

Signal test, using controlled reagents and plates:
ECL, positive control (tolerance range): 19 354 – 26 185

Baseline ECL (tolerance range): 16 - 56

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>	05 FEB 2019