



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD Read Buffer T (4X) Surfactant Free
Catalog #/Size: R92TD - Series
Lot Number: Y0060030
Expiration Date: 31 OCT 2021

Signal test, using controlled reagents and plates:
ECL, positive control (tolerance range): 48609 - 72914

Baseline ECL (tolerance range): ≤ 158

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>	08 JAN 2019