



CERTIFICATE OF ANALYSIS

COA - 07848

Product Description: Diluent 100

Catalog #/Size: R50AA - Series

Lot Number: Y0270243

Expiration Date: 31 JAN 2019

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

pH: 7.4 ± 0.05

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>X Talavera</i>	03 FEB 2017
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	07 FEB 2017