



CERTIFICATE OF ANALYSIS

COC - 00840-A

Product Description: MSD[®] Diluent 101

Catalog #/Size: R51AD-7 / 700 mL, R51AD-3 / 50 mL

Lot Number: M1010024

Expiration Date: 30 NOV 2018

Statement: This product was manufactured according to approved MSD documents.

Storage conditions: $\leq -10^{\circ}\text{C}$

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>Talavera</i>	19 OCT 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	19 OCT 2016