



CERTIFICATE OF CONFORMANCE

COC - 10484

Product Description: MSD[®] Diluent 2

Catalog #/Size: R51BB-2/200 mL; R51BB-3/40 mL; R51BB-4/8 mL

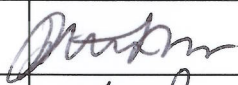
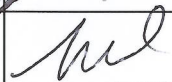
Lot Number: M0020077

Expiration Date: 31 AUG 2015

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	Meredith Yeager	Inventory Technician		01 May 2013
Review/Approval	Staci Leffel	Manufacturing Manager		01 MAY 2013