



CERTIFICATE OF CONFORMANCE

COC - 10450

Product Description: Diluent 101

Catalog #/Size: R51AD-7/ 700 mL

Lot Number: M1010006

Expiration Date: 31 OCT 2014

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		07/09/2012
Review/Approval	Staci Leffel	Manufacturing Manager		07/06/2012