



CERTIFICATE OF CONFORMANCE

COC - 10370

Product Description: MSD[®] Diluent 2

Catalog #/Size: R51BB-3/40 mL

Lot Number: M0020057

Expiration Date: 31 AUG 2013

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	Manufacturing Associate		09/14/2011
Review/Approval	Staci Leffel	Manufacturing Manager		09/08/2011
Review/Approval	Richard Mann	Quality Assurance		09/14/2011