



CERTIFICATE OF CONFORMANCE

COC - 10335

Product Description: MSD® Diluent 2
Catalog #/Size: R51BB-3/40 mL; R51BB-2/200 mL
Lot Number: M0020050
Expiration Date: 31 MAY 2013
Statement: This product was manufactured according to approved MSD documents.

Storage conditions: ≤ -10°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	<i>Meredith Yeager</i>	03/01/2011
Review/Approval	Staci Leffel	Manufacturing Manager	<i>Staci Leffel</i>	03/04/2011
Review/Approval	Richard Mann	Quality Assurance	<i>Richard M</i>	03/01/2011