



CERTIFICATE OF CONFORMANCE

COC - 10344

Product Description: MSD[®] Diluent 2
Catalog #/Size: R51BB-4/8 mL; R51BB-3/40 mL; R51BB-2/200 mL
Lot Number: M0020056
Expiration Date: 31 JUL 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		03/31/2011
Review/Approval	Staci Leffel	Manufacturing Manager		03/30/2011
Review/Approval	Richard Mann	Quality Assurance		03/31/2011