



CERTIFICATE OF CONFORMANCE

COC - 10259

Product Description: MSD® Diluent 2
Catalog #/Size: R56BB-3/50 mL; R51BB-2/200 mL
Lot Number: M0020041
Expiration Date: 31 AUG 2012
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	Janice Kim	Manufacturing Associate	Jai JK	05/10/2010
Review/Approval	Stacie Leffel	Kit Mfg Mng	Stacie Leffel	05/07/2010
Review/Approval	Richard Mann	Quality Assurance	Richard Mann	05/10/2010