



CERTIFICATE OF CONFORMANCE

COC - 10379

Product Description: MSD[®] Diluent 2

Catalog #/Size: R51BB-3/40 mL; R51BB-4/8 mL

Lot Number: M0020049

Expiration Date: 31 MAR 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	Rolanda Spratley	Quality Control		11 Oct 11
Review/Approval	Staci Leffel	Manufacturing Manager		10/11/2011
Review/Approval	Richard Mann	Quality Assurance		10/11/2011