



CERTIFICATE OF CONFORMANCE

COC - 10380

Product Description: MSD[®] Diluent 3

Catalog #/Size: R51BA-4/5 mL; R51BA-5/25 mL

Lot Number: M0030039

Expiration Date: 30 APR 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		11/21/11
Review/Approval	Staci Leffel	Manufacturing Manager		10/11/2011
Review/Approval	Richard Mann	Quality Assurance		12/11/2011