

CERTIFICATE OF COMPLIANCE

COC-10114

Product Description: MSD[®] Diluent 2

Catalog #/Product Size: R51BB-4/8 mL, R51BB-3/40 mL, or R51BB-2/200 mL

Lot Number: M0020033

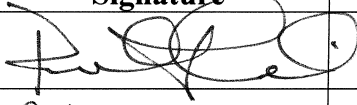

Expiration Date: March 31, 2012

Statement: This diluent was manufactured at MSD[®] according to MSD[®] SOPs and specifications.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

Prepared By:	Meagan McNamee	Date:	October 9, 2009
---------------------	----------------	--------------	-----------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Paul Goddwin		10/13/2009
Quality Assurance	Richard Mann		10/14/2009