

CERTIFICATE OF COMPLIANCE

COC-10113

Product Description: MSD[®] Diluent 3

Catalog #/Product Size: R51BA-4/5 mL, R51BA-5/25 mL, or R51BA-6/125 mL

Lot Number: M0030024

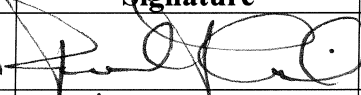

Expiration Date: 31 MAR 2012

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 5, 2009
---------------------	----------------	--------------	-----------------

Reviewed and Approved By

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