

CERTIFICATE OF COMPLIANCE

COC-10076

Product Description: MSD[®] Diluent 3

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

Lot Number: M0030014

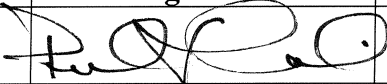

Expiration Date: February 2012

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:	September 8, 2009
---------------------	----------------	--------------	-------------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Paul Goodwin		9/10/09
Quality Assurance	Richard Mann		09/11/2009