

CERTIFICATE OF COMPLIANCE

COC-10118

Product Description: MSD[®] Diluent 2

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

Lot Number: M0020019

Expiration Date: February 2010

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 28, 2009
---------------------	----------------	--------------	--------------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Staci Laffel	<i>Staci Laffel</i>	10/04/2009
Quality Assurance	Richard Mann	<i>Richard Mann</i>	10/02/2009

REVISION HISTORY