

CERTIFICATE OF COMPLIANCE

COC-10188

Product Description: MSD[®] Diluent 2

Catalog #/Product Size: R51BB-2/200 mL

Lot Number: M0020038

Expiration Date: 30 APR 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:	November 25, 2009
---------------------	----------------	--------------	-------------------

Reviewed and Approved By

Function	Name (Print)	Signature	Date
Operations	Staci Leffel	<i>Staci Leffel</i>	11/25/2009
Quality Assurance	Richard Mann	<i>Richard Mann</i>	12/21/2009