

**CERTIFICATE OF COMPLIANCE**

**COC-10075**

**Product Description:** MSD<sup>®</sup> Diluent 2

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

**Lot Number:** M0020023



**Expiration Date:** February 2012

**Statement:** This diluent was manufactured at MSD according to MSD<sup>®</sup> SOPs and specifications.

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

<b>Prepared By:</b>	Meagan McNamee	<b>Date:</b>	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