

**CERTIFICATE OF COMPLIANCE**

**COC-10055**

**Product Description:** MSD<sup>®</sup> Diluent 3

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

**Lot Number:** M0030011

**Expiration Date:** February 2010

**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>	July 07, 2009
---------------------	----------------	--------------	---------------

**Reviewed and Approved By**

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

**REVISION HISTORY**