

## CERTIFICATE OF COMPLIANCE

COC-10049

**Product Description:** MSD<sup>®</sup> Human Serum Cytokine Assay Diluent

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

**Lot Number:** M0020016



**Expiration Date:** January 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>	<b>Meagan McNamee</b>	<b>Date:</b>	<b>June 10, 2009</b>
---------------------	-----------------------	--------------	----------------------

**Reviewed and Approved By**

Function	Name (Print)	Signature	Date
OPERATIONS	PAUL GOODWIN		6/10/2009
QA	Richard Mann		06/10/2009