

## CERTIFICATE OF COMPLIANCE

COC-10056

**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:** M0020020

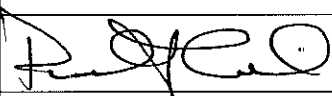

**Expiration Date:** February 2010

**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	PAUL GOODWIN		07/07/2009
Quality Assurance	Richard Mann		07/07/2009