

CERTIFICATE OF COMPLIANCE

COC-10050

Product Description: MSD[®] Human Serum Cytokine Antibody Diluent

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

Lot Number: M0030008



Expiration Date: December 2009

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:	June 10, 2009
---------------------	----------------	--------------	---------------

Reviewed and Approved By

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