





**CERTIFICATE OF CONFORMANCE**

COC - 00726

**Product Description:** MSD<sup>®</sup> Diluent 43  
**Catalog #/Size:** R50AG-2/50 mL; R50AG-1/10 mL  
**Lot Number:** M0430009  
**Expiration Date:** 30 NOV 2016  
**Statement:** This product was manufactured according to approved MSD documents.

Storage conditions: ≤-10°C

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

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Meredith Hutchinson	Inventory Technician		23 SEP 2014
<b>Review/Approval</b>	Latrice Corsey	Manufacturing Lead		23 SEP 2014