



---

**CERTIFICATE OF CONFORMANCE**

COC - 00783

**Product Description:** MSD® Diluent 43

**Catalog #/Size:** R50AG-2/50 mL; R50AG-1/10 mL



**Lot Number:** M0430012

**Expiration Date:** 30 JUN 2017

**Statement:** This product was manufactured according to approved MSD documents.

Storage conditions:  $\leq -10^{\circ}\text{C}$

**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		24 AUG 2015
<b>Review/Approval</b>	Staci Leffel	Manufacturing Manager		24 AUG 2015