





CERTIFICATE OF CONFORMANCE

COC - 00703

Product Description: MSD[®] Diluent 43
Catalog #/Size: R50AG-2/50 mL; R50AG-1/10 mL
Lot Number: M0430008
Expiration Date: 31 AUG 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.

	Name	Function	Signature	Date
Preparer	Meredith Hutchinson	Inventory Technician		07 MAY 2014
Review/Approval	Staci Leffel	Manufacturing Manager		08 MAY 2014