



CERTIFICATE OF ANALYSIS

COA - 04498

Product Description: MSD[®] Diluent 17

Catalog #/Size: R50KA-3/30 mL; R50KA-4/6 mL



Lot Number: M0170049

Expiration Date: 30 JUN 2016

Statement: This product was manufactured and tested for interferences 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.

	Name	Function	Signature	Date
Preparer	Meredith Hutchinson	Inventory Technician		04 JUN 2014
Review/Approval	Staci Leffel	Manufacturing Manager		06 JUN 2014