



CERTIFICATE OF ANALYSIS

COA - 06628

Product Description: MSD[®] Diluent 17

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



Lot Number: M0170052

Expiration Date: 31 OCT 2017

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	Production Control		07 FEB 2017
Review/Approval	Vanessa Castleberry	Quality Assurance		07 FEB 2017