



CERTIFICATE OF ANALYSIS

COA - 04713

Product Description: MSD Blocker A



Catalog #/Size: R93BA-2 / 250 mL

Lot Number: Y0180988

Expiration Date: 31 DEC 2016

Statement: This product was manufactured and tested for interferents according to approved MSD documents.

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		28 AUG 2014
Review/Approval	Anastasia Belilovets	Quality Control		28 AUG 2014