



## CERTIFICATE OF ANALYSIS

COA - 04420

**Product Description:** MSD Blocker A



**Catalog #/Size:** R93BA-2 / 250 mL

**Lot Number:** Y0180948

**Expiration Date:** 30 APR 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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		12 MAY 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		14May2014