



## CERTIFICATE OF ANALYSIS

COA - 11099

**Product Description:** MSD Blocker A



**Catalog #/Size:** R93BA-4 / 15 g

**Lot Number:** Y0180900

**Expiration Date:** 30 NOV 2014

**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		08 Nov 12
Review/Approval	Michael Eager	Quality Control		14 JAN 2013