



CERTIFICATE OF ANALYSIS

COA - 04415

Product Description: MSD Blocker A



Catalog #/Size: R93BA-4 / 15 g

Lot Number: Y0180961

Expiration Date: 31 AUG 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		09 MAY 2014
Review/Approval	Michael Eager	Quality Control		14May2014