



CERTIFICATE OF ANALYSIS



COA - 03596

Product Description: MSD[®] Mouse IL-13 Antibody
Catalog #/Size: D22AS-2 / 1 Plate; D22AS-3 / 5 Plates
Lot Number: D0032546
Expiration Date: 31 MAR 2014

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Source (host/species): Goat
Volume: 75 μ L per plate
Concentration: 50X
L/P: 7.41
Diluent: MSD Diluent 100
Azide added: 0.05% azide
Storage conditions: 2-8 $^{\circ}$ C, Dark

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		25 APR 2013
Review/Approval	Michael Eager	Quality Control		24 MAR 2013