



CERTIFICATE OF ANALYSIS

COA - 04096

Product Description: MSD[®] Human TNF- α Calibrator

Catalog #/Size: C01BH-2 / 1 Plate

Lot Number: A0031076

Expiration Date: 15 Months From Date of Receipt

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

Source (host/species): *E. coli*



Concentration: 1 μ g/mL

Purity: > 97% as determined by SDS-PAGE and visualized by silver stain.

Storage buffer: Diluent 1

Storage conditions: $\leq -70^{\circ}$ C

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 JAN 2014
Review/Approval	Michael Eager	Quality Control		10 FEB 2014