



## CERTIFICATE OF ANALYSIS

COA - 04097

**Product Description:** MSD<sup>®</sup> Human IL-6 Calibrator

**Catalog #/Size:** C01AK-2 / 1 Plate

**Lot Number:** A0031341

**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 µg/mL

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

Storage buffer: Diluent 1

Storage conditions: ≤ -70° 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