



CERTIFICATE OF ANALYSIS

COA - 03526

Product Description: MSD[®] Human M-CSF Calibrator

Catalog #/Size: C01IQ-2 / 1 Plate

Lot Number: A0031244

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): Recombinant protein derived from DNA sequence expressed in *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 to -20° 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		05 APR 2013
Review/Approval	Michael Eager	Quality Control		05 APR 2013