



CERTIFICATE OF ANALYSIS

COA - 04197

Product Description: MSD[®] Human MMP 3-Plex Calibrator Blend

Catalog #/Size: C0034-2 / 1 Plate

Lot Number: A0031457

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): MMP-1, MMP-3: NS0; MMP-9: CHO



Concentration: MMP-1, MMP-3: 1µg/mL; MMP-9: 5µg/mL

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

Storage buffer: TCNB Additive

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 FEB 2014
Review/Approval	Michael Eager	Quality Control		04 MAR 2014