



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0028

**Expiration Date:** 31 DEC 2021

**Storage Condition:** 2-8°C

Analyte Description	Amount of Analyte in Calibrator (pg)	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) (pg/mL)	Concentration of Standard 1 (after 5-fold Dilution) (pg/mL)
Hu Eotaxin-2	8 375	33 500	6 700
Hu GRO-α	3 038	12 150	2 430
Hu I-309	3 450	13 800	2 760
Hu MCP-2	2 738	10 950	2 190
Hu MCP-3	5 500	22 000	4 400
Hu M-CSF	2 513	10 050	2 010
Hu MIF	37 500	150 000	30 000
Hu MIP-5	37 875	151 500	30 300
Hu TRAIL	13 875	55 500	11 100
Hu YKL-40	6 225	24 900	4 980

**Statement:** The above calibrator lot was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	22 FEB 2017