



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Proinflammatory Panel 1 (rat) Control Pack

**Catalog Number:** C4044-1

**Lot Number:** CP000074

**Expiration Date:** 30 APR 2023

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

Units = pg/mL	Assay								
	rt IFN- $\gamma$	rt IL-1 $\beta$	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- $\alpha$
Reconstituted Control 1 (A00C0582)	5130	12 000	977	3640	10 600	1180	23 300	1370	868
Reconstituted Control 2 (A00C0583)	1550	3180	275	1820	3480	569	7020	400	266
Reconstituted Control 3 (A00C0584)	454	1090	84.9	927	1130	289	2120	123	87.7

**Reconstitution Volume:** 250  $\mu$ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0582	2-8°C	30 APR 2023
Proinflammatory Panel 1 (rat) Control 2	A00C0583	2-8°C	30 APR 2023
Proinflammatory Panel 1 (rat) Control 3	A00C0584	2-8°C	30 APR 2023

Statement: This product 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>	21 JUN 2018