

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Acute Phase Protein Panel 1 (rat) Kit
Kit Catalog Numbers: K15175C-series
Kit Lot Number: K0040425
Expiration Date: 30 Apr 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	AGP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	A2M	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	AGP	0.150	80% - 120%	< 20%	150	75% - 125%	< 25%
4	A2M	1.00	80% - 120%	< 20%	400	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	AGP	< 1.0%
4	A2M	< 1.0%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	AGP	20X Calibrator	18700	ng/mL
		Control 1	234	ng/mL
		Control 2	58.4	ng/mL
		Control 3	14.6	ng/mL
4	A2M	20X Calibrator	12800	ng/mL
		Control 1	160	ng/mL
		Control 2	40	ng/mL
		Control 3	10	ng/mL

Control data was generated utilizing internal reference controls.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

CERTIFICATE OF ANALYSIS

Page 2 of 2

Standard Curve Data:


AGP - Spot 1			A2M - Spot 4		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
935	97,600	2.1	640	277,000	1.8
234	75,200	2.1	160	121,000	3.3
58.4	42,100	3.4	40.0	33,700	2.8
14.6	15,700	1.4	10.0	8,220	2.2
3.65	4,390	1.2	2.50	2,210	2.6
0.913	1,230	2.1	0.625	698	2.6
0.228	417	2.7	0.156	340	2.6
0.00	152	8.5	0.00	214	5.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Acute Phase Protein Panel 1 (rat) Plate	Z0056193	2-8°C	31 Aug 2020
Anti-rat AGP Antibody	D0040763	2-8°C	31 Aug 2021
Anti-rat A2M Antibody	D0040762	2-8°C	31 Aug 2021
Acute Phase Protein Panel 1 (rat) Calibrator Blend	A0040205	≤ -70°C	30 Apr 2020
Diluent 100	Y0270253	2-8°C	30 Apr 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		06 MAR 2019