



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Acute Phase Protein Panel 1 (rat) Calibrator Blend

Catalog Number: C0175-2

Lot Number: A0040219

Expiration Date: 31 MAR 2021

Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration of Materials Tested:

	Assay	
	AGP	A2M
Calibrator (ng/mL)	8760	5540

Concentration: 20X

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	15 NOV 2018