



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07935

**Product Description:** Rat BNP Controls

**Catalog Number:** C01XX-2

**Lot Number:** CP000046

**Expiration Date:** 28 FEB 2022

### Concentration of Materials Tested:

Units = pg/mL	Rat BNP
Description	
Rat BNP Control 1	924
Rat BNP Control 2	204
Rat BNP Control 3	52.2

Rat BNP Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Rat BNP Control 1	A00C0457	$\leq -70^{\circ}\text{C}$	28 FEB 2022
Rat BNP Control 2	A00C0458	$\leq -70^{\circ}\text{C}$	28 FEB 2022
Rat BNP Control 3	A00C0459	$\leq -70^{\circ}\text{C}$	28 FEB 2022

**Fold Dilution:** 2X

**The above product is intended for Internal Testing Use Only. Not for use by Customers.**

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	06 MAR 2017