



Certificate of Analysis / Data Sheet

Product: Vancomycin Hydrochloride**Catalog No: RC-206****Release Date: 8/25/20****Size: 5g****Retest Date: 8/25/24****Lot No: 202508****Storage: Desiccated at -20°C****Specifications****CAS Number: 1404-93-9****Appearance: White to off-white or tan to brown**

Test	Specification	Result
Identification (IR)	Conforms	PASS
Identification (Chloride)	Conforms	PASS
Solubility	Soluble in water, insoluble in ether and chloroform	PASS
pH (50mg/ml water)	2.5-4.5	3.11
Water	≤ 5.0%	2.95%
Residue on Ignition	≤ 1.0%	0.03%
Assay (anhydrous basis)	≥ 900µg/mg	1028µg/mg
Vancomycin B	≥90.0%	96.1%
Related Substance:		
N-demethyl Vancomycin B	≤ 4.0%	0.83%
Desamido Vancomycin B	≤ 4.0%	0.08%
Agluco Vancomycin B	≤ 4.0%	0.05%
Desvancosaminy Vancomycin B	≤ 4.0%	0.19%
Other	≤ 5.0%	0.67%
Total impurities	≤ 10.0%	3.85%
Monodechlorovancomycin	≤ 4.7%	1.3%



Bacterial Endotoxins	≤ 0.33 EU/mg	< 0.25 EU/mg
Total Aerobic Microbial Count	≤ 100 CFU/g	< 100 CFU/g
Combined Yeast & Mold Count	≤ 10 CFU/g	< 10 CFU/g

This document has been produced electronically and is valid without signature

Date of Print: September 14, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP