



Certificate of Analysis / Data Sheet

Product: X-Gluc (5-Bromo-4chloro-3indoxyl- β -D-glucuronic acid, CHX salt)
Catalog No: RC-215 **Release Date:** 9/2/20
Size: 1g **Retest Date:** 9/2/24
Lot No: 200209 **Storage:** Desiccated at -20°C

Specifications

CAS Number: 114162-64-0

Appearance: White to off-white crystalline powder

Test	Specification	Result
Color (1% in DMF/water 1:1)	Colorless	Colorless
Turbidity (1% in DMF/water 1:1)	≤ 4.0 NTU	0.3 NTU
Water Content (K. Fischer)	2.5 – 4.5%	3.4%
Rotation	-91.0 to -87.0	-88.9
Perchloric Acid Titration (0.1M)	99.0 – 101.0%	99.8%
Purity (HPLC)	≥ 99.0 area-%	99.4 area-%
TLC	Corresponds	PASS
FT-IR Spectrum	Corresponds	PASS

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

