



Certificate of Analysis

Product: Octyl -beta-D-Glucopyranoside

Catalog #: DG015

Size: 1g

Lot#: 111909

Molecular Formula: C₁₄H₂₈O₆

Molecular Weight: 292.38

Tests:

Appearance: White crystals	Passed	Yes / No
Purity (Min 98.0%)	Passed	Yes / No
Specific Rotation (Min -32.0 degrees)	Passed	Yes / No
Protease (None)	Passed	Yes / No
Solubility (10%, water)	Passed	Yes / No

Storage: Store at -20⁰C and keep dry. Warm it to room temperature before opening.

This document has been produced electronically and is valid without signature

Date of Print: December 8, 2011

G-Biosciences/ Genotech

92 Weldon Parkway, St. Louis, MO 63043, USA
Ph: 314-991-6034, Fax: 314-991-1504

Octyl beta Glucopyranoside DG015 C of A/Com#3/Rev 04.19.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com