



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: PEG 8000

Catalog No: RC-078

Release Date: 7/30/20

Size: 1 kg

Retest Date: 7/30/26

Lot No: 203007

Storage: Room Temp.

Specifications

CAS Number: 25322-68-3

Appearance: Fine white powder

Test	Specification	Result
Hydroxyl number	13.6	PASS
Melting point	58-59°C	PASS

This document has been produced electronically and is valid without signature

Date of Print: September 2, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com