



Certificate of Analysis

Product: 1,5-Diphenylcarbazine

Catalog #: BTNM-0028

Quantity: 25g

Lot#: 181603

CAS number: 140-22-7

MW: 242.28

Appearance: Pale grey powder

Tests:

	Actual	Disposition
Purity (HPLC)	99%	Passed
Melting Point (Degree C)	171.1-172.2	Passed
Identification (FTIR)	Conforms	Passed

Storage Temp: Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

Date of Print: April 17, 2018

G-Biosciences / Genotech

9800 Page Avenue, St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

