



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Nuclear Fast Red Solution (Kernechtrot)

Catalog No: 786-1053

Release Date: 01/04/21

Size: 100mL

Retest Date: 01/04/25

Lot No: 210401

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, red liquid

Test	Specification	Result
Stain	Pink to rose color	PASS
O.D. @ 535nm	0.560-0.850	0.845

This document has been produced electronically and is valid without signature

Date of Print: February 23, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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