



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: 786-1055

Release Date: 8/17/20

Size: 1 L

Retest Date: 8/17/24

Lot No: 201708

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	0.776

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