



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Cytoscan LDH Cytotoxicity Assay kit

Catalog No: 786-324

Release Date: 04/28/2023

Size: Kit

Retest Date: 04/23/2029

Lot No: 232504

Storage: store as recommended

Specifications

CAS Number: not applicable

Appearance: not applicable

Test

Specification

Result

This kit performs as per corresponding
and/or companion protocol

Yes/No

PASS

This document has been produced electronically and is valid without signature

Date of Print: June 9, 2023

G-Biosciences/Geno Technology, Inc.

Rev 02/13/2022 DB



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