



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** Cytoscan LDH Cytotoxicity Assay kit

**Catalog #:** 786-324

**Size:** Kit

**Lot#:** 202901

**Re-test date:** 01/2025

### Analysis

#### Passes

This kit performs as per corresponding and/or companion protocol

Yes/~~No~~

This document has been produced electronically and is valid without signature

**Date of Print:** March 24, 2020

*For Research Use Only*

### **G-Biosciences / Genotech**

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)