



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

**Catalog #:** 786-324

**Size:** Kit

**Lot#:** 210212

**Re-test date:** 12/2031

### 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:** May 4, 2022

*For Research Use Only*

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

9800 Page Avenue

St. Louis, MO 63132, USA

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

Aprotinin C of A/Rev 04.29.15-IA



think proteins! think G-Biosciences!

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