



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: Bacterial Protease Arrest

Cat #: 786-330

Lot #: 211106

Re-Test date: 06/2026

TESTS:

Tests performed using Protease Arrest, Cat # 786-108. protocol. Protease inhibition is performed using G-Biosciences's protease preparation; pancreas extract. Protease inhibition is analyzed by measuring activity of dye-labeled protein as substrate. Incubation time 3 hours. Triplicate of each sample analyzed.

Protease Arrest at 1X concentration

Control (average O.D.)	0.388
Test, Inhibition (Average O.D.)	0.013
% Inhibition	96.6%

Specification for Pass: For the assay to pass, the % of inhibition must be > 90%

This document has been produced electronically and is valid without a signature

Date of print: January 20, 2022

G-Biosciences / Genotech

9800 Page Avenue, St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Protease Arrest CA/Com#3, Rev 12.22.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com