



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

Certificate of Analysis

Product: Phosphatase Arrest-I

Cat #: 786-450

Lot #: 211011

Re-test date: 11/2026

TESTS:

Tests performed using Phosphatase Arrest, Cat # 786-450 protocol. Phosphatase inhibition was performed using *G-Biosciences's* phosphatase preparation and was analyzed by measuring the absorbance of control and test samples at 400nm.

Phosphatase Arrest-I at 1X concentration:

Control (average OD)	0.992
Test Sample (average O.D.)	0.275
% Inhibition	78.7%

Specification for Pass: For the assay to pass, the % of inhibition must be > 70%

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

Date of print: December 21, 2021

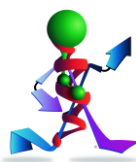
G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

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

Phosphatase Arrest-I C of A/Com#3 Rev 01.16.12/IA



think proteins! think G-Biosciences!

www.GBiosciences.com