



Certificate of Analysis / Data Sheet

Product: Phosphatase Arrest III

Catalog No: 786-452

Release Date: 12/12/19

Size: 1ml

Retest Date: 12/12/24

Lot No: 191212

Storage: 4°C

Specifications

Appearance: Clear solution

Tests performed using Phosphatase Arrest-III, Cat # 786-452 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. The following tests were performed with Phosphatase Arrest III at 1X concentration.

Test	Specification	Result
Control (avg. OD)	N/A	0.484
Test Sample (avg. OD)	N/A	0.047
% Inhibition	> 80%	90.3%

This document has been produced electronically and is valid without signature

Date of Print: December 24, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

