



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



Certificate of Analysis

Product: SAM510: SAM Methyltransferase Assay

KIT COMPONENTS

Cat # 786-430

SAM Methyltransferase Assay Buffer	20ml
SAM Methyltransferase Assay Buffer Additive	0.2ml
SAM510 Enzyme Mix	5 x 300µl
SAM Colorimetric Mix	1
Positive Control: Adenosylhomocysteine 1mM]	0.2ml
S-Adenosylmethionine	5 x 200µl
HCl Assay Reagent	1 ml
Dry Ice	1

Lot#: 192301

Tests:

SAM Methyltransferase Assay

Passed

Yes / ~~No~~

SAM510 Assay quantitatively assays the Positive Control: Adenosylhomocysteine (AdoHcy). 0-100µM AdoHcy was assayed with SAM510.

Storage Conditions: -70°C

This document has been produced electronically and is valid without signature

Date of Print: January 14, 2022

G-Biosciences / Genotech

9800 Page Ave

St. Louis, MO 63132, USA

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

SAM510: SAM Methyltransferase Assay C of A/02.05.09-IA/Common



think proteins! think G-Biosciences!

www.GBiosciences.com