



Certificate of Analysis / Data Sheet

Product: FCCP [Carbonyl cyanide 4-(trifluoromethoxy) phenylhydrazone]
Catalog No: 786-1317 **Release Date:** 2/12/19
Size: 50mg **Retest Date:** 6/15/19
Lot No: 192501 **Storage:** -20°C

Specifications

CAS Number: 370-86-5

Appearance: Yellow powder

Test	Specification	Result
HPLC	≥ 98%	100.0%
Mass spec	Report result	M-H ⁺ : 253.1
TLC	≥ 98%	100%
UV	Report result	λ max: 355, 239 nm

This document has been produced electronically and is valid without signature

Date of Print: December 20, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

