



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: PMSF

Catalog No: 786-055

Size: 5g

Lot No: 181105A

Release Date: 08/29/2018

Retest Date: 7/20/2019

Storage: RT

Specifications

CAS Number: 329-98-6

Appearance: White Needle-like Crystals

Test	Specification	Result
Loss on Drying	≤ 0.5%	< 0.5%
Melting Point	94.0-94.0°C	93.1-93.9°C
Purity (GC)	≥ 98.5%	99.67%
Solubility	Clear, Colorless Solution in Chloroform and Diethyl Ether	Passed
TLC Spot	One Spot	Conforms

This document has been produced electronically and is valid without signature

Date of Print: January 15, 2019

G-Biosciences/Geno Technology, Inc.
9800 Page Avenue, St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Rev 02/13/17 CCH



think proteins! think G-Biosciences www.GBiosciences.com