



Certificate of Analysis

AMP+ Mass Spectrometry Kit

Item #: 710000
Batch #: 0545448

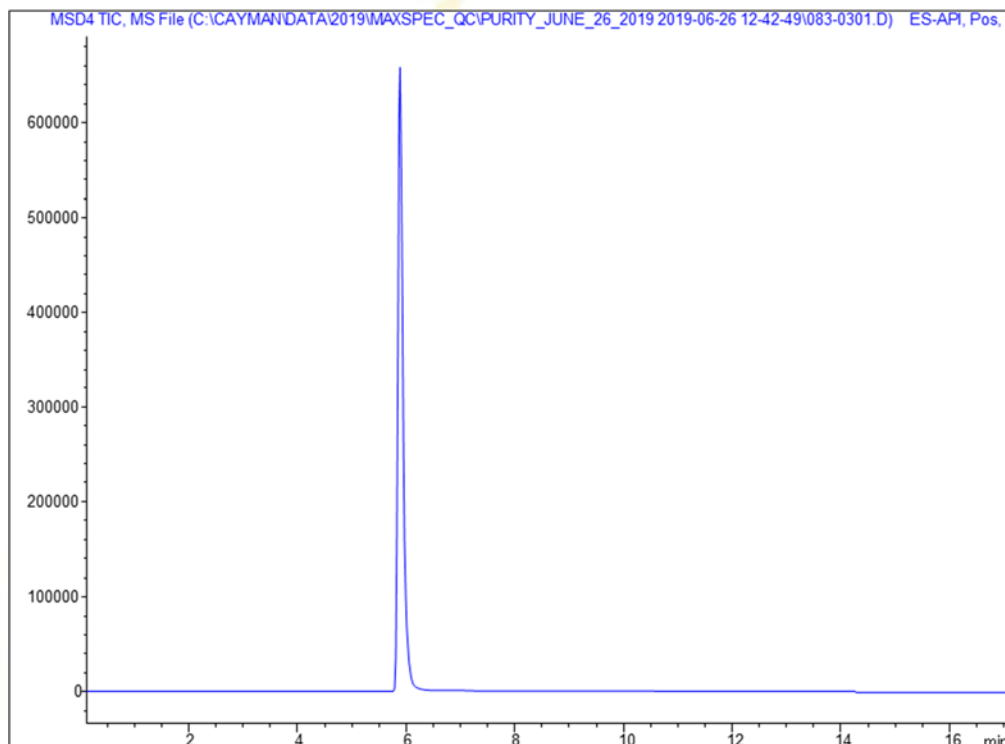
Contents

Component	Item / Batch #	Description
AMP+ Derivatizing Reagent	710001 / 0487912	20mM solution in acetonitrile
HOBt (1-Hydroxybenzotriazole monohydrate)	710002 / 0487913	20mM solution in 99:1 acetonitrile:DMF
EDC (1-Ethyl-(3-dimethylaminopropyl)carbodiimide hydrochloride)	710003 / 0487914	640mM aqueous solution
Acetonitrile / DMF solution	710004 / 0487915	4:1 acetonitrile:DMF solution
Autosampler Vial Inserts	710005 / 0487916	200uL plastic polyspring inserts

Quality Information

Qualifier	Method	Limit	Result	Meets Specification
AMP Derivatizing Reagent	NMR	Conforms	Conforms	Y
HOBt	Purity (assay)	≥99%	100%	Y
EDC	Purity (assay)	≥99%	99.9%	Y

AMP+ Derivatized PGF_{2α}



Agilent 1100 HPLC / Agilent 6120
Single Quadrupole MS

Sample: 8μg PGF_{2α} derivatized
with AMP+ according to kit
specifications

Method:

A: H₂O + 0.1% Formic Acid

B: MeCN + 0.1% Formic Acid

0min: 10%B

1min: 10%B

8min: 98%B

11min: 98%B

11.1min: 10%B

17min: 10%B

Phenomenex Kinetex EVO C18,
3.0x50mm, 2.6μ,
400μl/min, 30°C

Cayman Chemical Co.

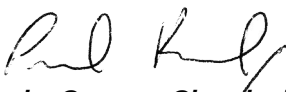
1180 E. Ellsworth Rd
Ann Arbor, MI 48108
www.caymanchem.com

Contact Us:

1-800-364-9897 or (734)971-3335
Fax: (734)971-3640
Email: techserv@caymanchem.com

For R&D purposes only. Not intended for human or animal use.

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiration date when stored unopened as recommended.

Approval: 
Paul Kennedy, Cayman Chemical Director of Analytical Chemistry

Certification Date: 16NOV2018

Revision History

Revision #	Date	Reason for Revision
01	16NOV2018	Initial version

Disclaimers**Material Safety Data**

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some but not all of the information required for the safe and proper use of this material. Before use, review the complete Material Safety Data Sheet, which has been sent via email to your institution.

Warranty and Limitation of Remedy

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specification.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver of Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitations of Remedy located on our website and in our catalog.

This Certificate shall not be reproduced except in full, without written approval from the Cayman Chemical Director of Analytical Chemistry.

