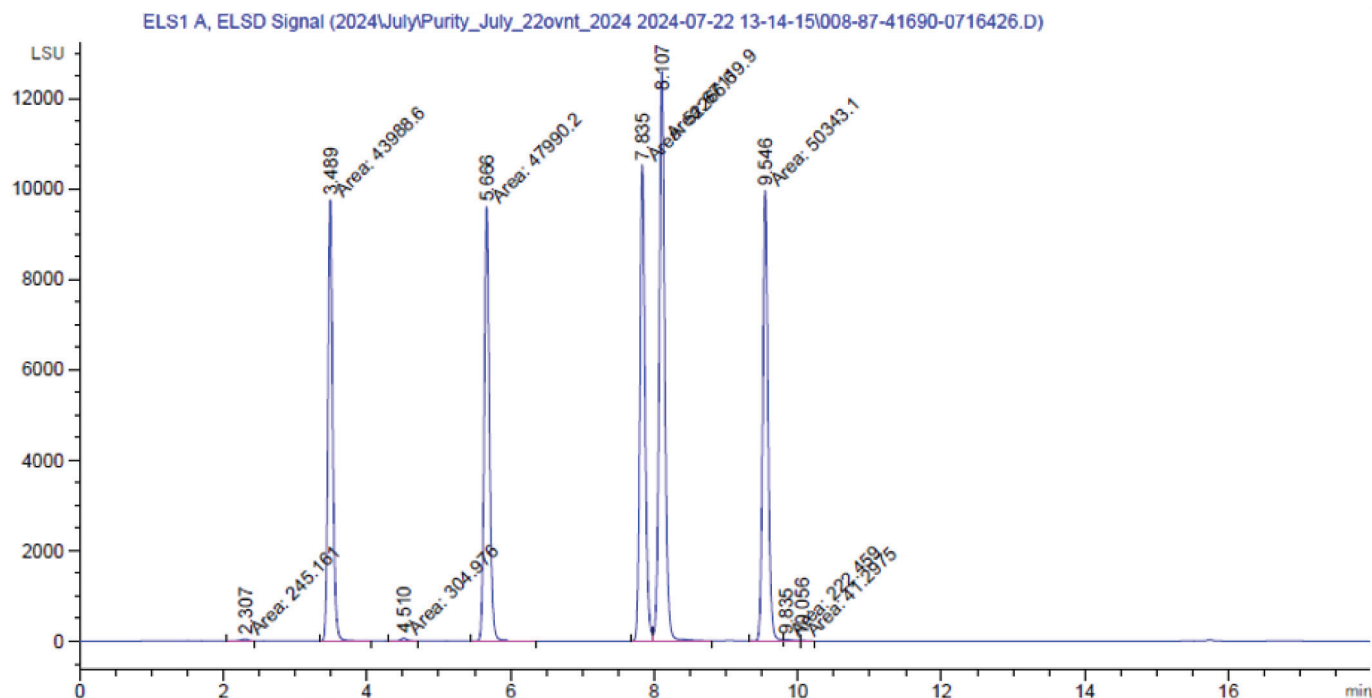


Quality Information

Test	Specification	Result
HPLC-ELSD	Identities confirmed	Pass



E. Johnson

Certification Date: 31JUL2024

Page 1 of 3

CERTIFICATE of ANALYSIS



Quality Control Table

Item No.	Item Name	RT (min)
27281	1,2,3-Trihexanoyl Glycerol	3.5
26952	1,2,3-Trioctanoyl Glycerol	5.7
26958	1,2,3-Tridecanoyl Glycerol	7.8
23336	1,2,3-Trilauroyl Glycerol	9.5
10030659	(+)- α -Tocopherol acetate	8.1
HPLC Conditions: Agilent 1260 Series HPLC-ELSD		
Mobile Phase A: 60:40:0.1:10 mM Water:Acetonitrile:Formic Acid:Ammonium Formate		
Mobile Phase B: 90:10:0.1:10 mM Water:Acetonitrile:Formic Acid:Ammonium Formate		
Column: Gemini C18 5 μ m 4.6 x 150 mm		Flow Rate: 1 ml/min
Gradient: 0-10 min 50-90% B; 10-13 min hold 99% B; 13-13.1 min 99%B to 50%B; 13.1-18 min ReEQ 50% B		

Mixture was prepared using gravimetric solution preparation procedures from neat analytical standards. The concentration of this mixture was prepared with an estimated error of $\pm 10\%$.

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	31JUL2024	Initial version
02	22AUG2024	Updated HPLC conditions

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 Company.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com