

CERTIFICATE of ANALYSIS



MaxSpec[®] Reference Plasma (human)

Item No.:	42573
Batch No.:	0802194
Expiry Date:	14MAR2030 (valid from date of certification)
Contents:	1 ml (overfilled to 1.1 ml) pooled human plasma (in lithium heparin), filtered (0.2 µm), and provided in 2 ml screw cap vials.
Storage & Handling:	Store upright and unopened at -80°C. <i>Biohazard statement: Please use universal precautions when handling this material. Material should be handled as if capable of transmitting infectious agents. No test method can provide total assurance that HIV, Hepatitis B, Hepatitis C, or other infectious agents are absent. All biological products should be handled at the Biosafety Level 2 as recommended by the CDC/NIH manual "Biosafety in microbiological and biomedical laboratories, from potentially infectious human serum or blood specimens." This product is being sold for research purposes only and is not intended as a source of material for clinical or diagnostic use.</i>

Quality Control

Donors used for this product have been tested with FDA CBER licensed screening tests (Title 21 - CFR part 610.40) to show the following: Hepatitis B surface antigen negative, HBV NAT negative, HIV 1&2 antibody negative, HIV NAT negative, HCV antibody negative, HCV NAT negative, syphilis negative, and West Nile virus NAT negative. Donors have tested negative by a licensed test for antibodies to *Trypanosoma cruzi* either on the current donation or at least one previous donation.

Donor Information

Gender	Age
Male	26
Male	50
Male	61
Male	39
Male	36
Male	47
Male	22
Male	33
Female	32
Female	32
Female	62
Female	42
Female	28
Female	59
Female	26

Approval: 

Title: Manager, Analytical Chemistry

Certification Date: 14MAR2025

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 expiry date when stored unopened as recommended.

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	14MAR2025	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 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