

PRODUCT INFORMATION

MaxSpec[®] Reference Plasma (Human)

Item No. 42573

Supplied as: A 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: -80°C

Stability: ≥5 years; *Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.*

Special Conditions: Store upright and unopened at -80°C.

Description

MaxSpec[®] Reference Plasma (Human) contains 1 ml of pooled human plasma intended for use as a reference material for the analysis of lipids or metabolites. This product includes a batch-specific downloadable table of analytes with their associated quantitative values and references to applicable methods. The table will continue to be updated as new analytes are tested and added. This product has undergone significant validation and traceability testing and can be used as a quality control in lipidomics method development, performance validation, or as an interlab control material. For more information, please see Cayman's application note: Characterization of Human Plasma as a Reference Material for Lipid Analysis Using LC-MS.

The human plasma used for the isolation of this product was found to be negative for HIV 1/2 antibody, HIV (NAT), HBsAg, HBV (NAT), HCV antibody, HCV (NAT), syphilis, and West Nile virus (NAT) by FDA-licensed screening assays. In addition, the donor has tested negative via a licensed test for T. cruzi antibodies either on the current donation or at least one previous donation. Human blood products should always be properly handled in accordance with safe handling procedures. Disposal should follow state and federal regulations and is the responsibility of the customer.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/11/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

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

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM