

PRODUCT INFORMATION

Polyamines Supplement Mix (1,000X)

Item No. 43829

Synonym: Putrescine/Spermidine/Spermine Mix

Supplied as: A solution

Storage: -20°C

Stability: ≥2 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Polyamines Supplement Mix (1,000X) is a cell culture media supplement and a mixture of essential polyamine cations known to support cell growth. It is used as a component of the CEPT small molecule cocktail to enhance the viability of human pluripotent stem cells (hPSCs).^{1,2} Polyamines Supplement Mix (1,000X) is provided as a 1,000X concentrated stock solution in water and may be added directly to cell culture media. This product is also available as part of Cayman's Stem Cell Survival Reagent Kit (Item No. 502884).

References

1. Chen, Y., Tristan, C.A., Chen, L., *et al.* A versatile polypharmacology platform promotes cytoprotection and viability of human pluripotent and differentiated cells. *Nat. Methods* **18**(5), 528-541 (2021).
2. Ryu, S., Weber, C., Chu, P.-H., *et al.* Stress-free cell aggregation by using the CEPT cocktail enhances embryoid body and organoid fitness. *Biofabrication* **16**, 015016 (2024).

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, 09/12/2025

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