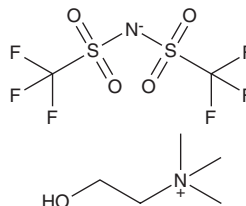


PRODUCT INFORMATION

Choline bis(trifluoromethylsulfonyl)imide

Item No. 43921

CAS Registry No.: 827027-25-8
Formal Name: 2-hydroxy-N,N,N-trimethyl-ethanaminium salt with 1,1,1-trifluoro-N-[(trifluoromethyl)sulfonyl]methanesulfonamide
Synonym: [Choline][NTf₂]
MF: C₅H₁₄NO • C₂F₆NO₄S₂
FW: 384.3
Purity: ≥98%
Supplied as: A low melting solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Choline bis(trifluoromethylsulfonyl)imide is supplied as a low melting solid. A stock solution may be made by dissolving the choline bis(trifluoromethylsulfonyl)imide in the solvent of choice, which should be purged with an inert gas. Choline bis(trifluoromethylsulfonyl)imide is soluble (≥10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of choline bis(trifluoromethylsulfonyl)imide can be prepared by directly dissolving the low melting solid in aqueous buffers. Choline bis(trifluoromethylsulfonyl)imide is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Choline bis(trifluoromethylsulfonyl)imide is an ionic liquid.¹ It has been used to extract phenolic compounds from an aqueous environment and as a thermally responsive ionic liquid in a closed-loop osmotic heat engine.^{1,2}

References

1. Cesari, L., Canabady-Rochelle, L., and Mutelet, F. Extraction of phenolic compounds from aqueous solution using choline bis(trifluoromethylsulfonyl)imide. *Fluid Phase Equilib.* **446**, 28-35 (2017).
2. Zhong, Y., Wang, X., Feng, X., et al. Osmotic heat engine using thermally responsive ionic liquids. *Environ. Sci. Technol.* **51(16)**, 9403-9409 (2017).

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, 10/06/2025

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