

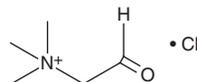
PRODUCT INFORMATION



Betaine Aldehyde (chloride)

Item No. 17270

CAS Registry No.: 7758-31-8
Formal Name: N,N,N-trimethyl-2-oxo-ethanaminium, monochloride
MF: C₅H₁₂NO • Cl
FW: 137.6
Purity: ≥95%
Supplied as: A crystalline 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

Betaine aldehyde (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the betaine aldehyde (chloride) in the solvent of choice, which should be purged with an inert gas. Betaine aldehyde (chloride) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

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 betaine aldehyde (chloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of betaine aldehyde (chloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Betaine aldehyde is the physiological intermediate in the oxidation of choline to betaine. This step is involved in glycine, serine, and threonine metabolism.

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, 11/17/2022

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