

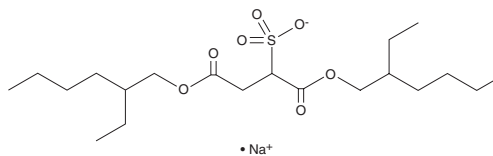
PRODUCT INFORMATION



Docusate (sodium salt)

Item No. 30692

CAS Registry No.: 577-11-7
Formal Name: 2-sulfo-butanedioic acid, 1,4-bis(2-ethylhexyl) ester, monosodium salt
Synonyms: Aerosol OT, AOT, Dioctyl sulfosuccinate, DOSS, KM 10-1610
MF: C₂₀H₃₇O₇S • Na
FW: 444.6
Purity: ≥95%
Supplied as: A 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

Docusate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the docusate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Docusate (sodium salt) is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of docusate (sodium salt) in these solvents is approximately 15 and 10 mg/ml, respectively.

Docusate (sodium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, docusate (sodium salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Docusate (sodium salt) has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Docusate is a surfactant.¹ It forms reverse micelles in oils and increases the solubility of water in oil by forming microemulsions. Docusate has been used in protein extraction, cryo-enzymology, the preparation of ceramic nanoparticles and gelatin-based nanopallets, and as a charge control agent in particle migration imaging.

Reference

1. De, T.K. and Maitra, A. Solution behaviour of aerosol OT in non-polar solvents. *Adv. Colloid Interfac.* **59**, 95-193 (1995).

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, 05/30/2024

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