

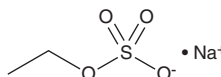
PRODUCT INFORMATION



Ethyl sulfate (sodium salt)

Item No. 39408

CAS Registry No.: 546-74-7
Formal Name: Sulfuric acid, monoethyl ester, monosodium salt
Synonyms: Ethyl sodium sulfate, Sodium ethyl sulfate
MF: $C_2H_5O_4S \cdot Na$
FW: 148.1
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Description

Ethyl sulfate (sodium salt) (Item No. 39408) is an analytical reference standard categorized as a metabolite of ethanol.¹ The presence of ethyl sulfate in urine or blood has been used as a biomarker for recent alcohol consumption. This product is intended for research and forensic applications.

Reference

1. Helander, A. and Beck, O. Ethyl sulfate: A metabolite of ethanol in humans and a potential biomarker of acute alcohol intake. *J. Anal. Toxicol.* **29(5)**, 270-274 (2005).

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, 03/03/2026

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