

PRODUCT INFORMATION



CENTA

Item No. 31534

CAS Registry No.: 80072-86-2
Formal Name: (6R,7R)-3-[[[3-carboxy-4-nitrophenyl]thio]methyl]-8-oxo-7-[[2-(2-thienyl)acetyl]amino]-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid

MF: C₂₁H₁₇N₃O₈S₃

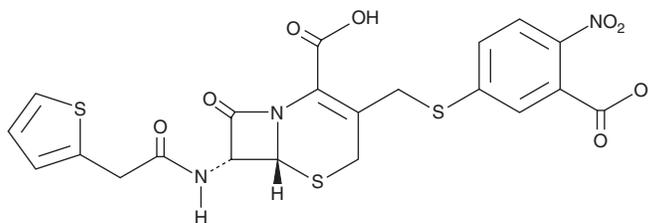
FW: 535.6

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

CENTA is supplied as a solid. A stock solution may be made by dissolving the CENTA in the solvent of choice, which should be purged with an inert gas. CENTA is soluble in the organic solvent methanol.

Description

CENTA is a colorimetric cephalosporin substrate for β -lactamases.^{1,2} Upon hydrolysis by β -lactamases, CENTA turns from light yellow to chrome yellow, which can be quantified by colorimetric detection at 405 nm as a measure of β -lactamase activity.

References

1. Jones, R.N., Wilson, H.W., Novick Jr., W.J., *et al.* In vitro evaluation of CENTA, a new beta-lactamase-susceptible chromogenic cephalosporin reagent. *J. Clin. Microbiol.* **15**(5), 954-958 (1982).
2. Bebrone, C., Moali, C., Rival, S., *et al.* CENTA as a chromogenic substrate for studying beta-lactamases. *Antimicrob. Agents Chemother.* **45**(6), 1868-1871 (2001).

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, 11/10/2022

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