

PRODUCT INFORMATION



Orlistat Degradation Product (sodium salt)

Item No. 27179

Formal Name: [2S-(2R*,3R*,5R*)]-2-hexyl-3,5-dihydroxy-hexadecanoic acid, monosodium salt

Synonym: Orlistat M3 Impurity

MF: C₂₂H₄₃O₄ • Na

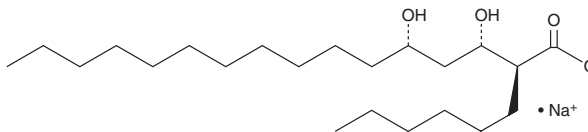
FW: 394.6

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

Orlistat degradation product (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the orlistat degradation product (sodium salt) in the solvent of choice, which should be purged with an inert gas. Orlistat degradation product (sodium salt) is soluble in the organic solvent methanol.

Description

Orlistat degradation product is a degradation product of the digestive lipase inhibitor orlistat (Item No. 10005426).¹ It is formed *via* hydrolytic and thermal degradation as well as digestion by human carboxyl ester lipase.^{1,2}

References

1. Stalder, H., Schneider, P.R., and Oesterhelt, D. Tetrahydrolipstatin: Thermal and hydrolytic degradation. *Helv. Chim. Acta.* **73(4)**, 1022-1036 (1990).
2. Stalder, H., Oesterhelt, G., and Borgström, B. Tetrahydrolipstatin: Degradation products produced by human carboxyl-ester lipase. *Helv. Chim. Acta.* **75(5)**, 1593-1603 (1992).

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, 12/21/2022

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