

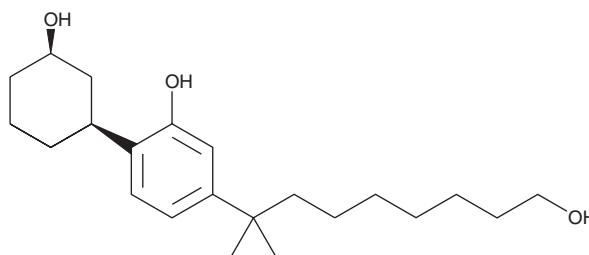
PRODUCT INFORMATION



CP 47,497-C8-homolog C-8-hydroxy metabolite

Item No. 9000773

CAS Registry No.: 1554485-48-1
Formal Name: *rel*-3-hydroxy-4-[(1*R*,3*S*)-3-hydroxycyclohexyl]- θ,θ -dimethyl-benzeneoctanol
MF: $C_{22}H_{36}O_3$
FW: 348.5
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 204, 218, 276 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥ 4 years



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

Description

CP 47,497-C8-homolog is similar to CP 47,497, a monophenol synthetic cannabinoid (CB) that potently imitates natural CBs. CP 47,497-C8-homolog avidly binds the CB_1 receptor ($K_i = 0.83$ nM).¹ It has been identified as an adulterant of herbal blends.² CP 47,497-C8-homolog C-8-hydroxy metabolite (Item No. 9000773) is a potential metabolite of CP 47,497-C8-homolog. This product is intended for research and forensic purposes.

References

1. Melvin, L.S., Milne, G.M., Johnson, M.R., *et al.* Structure-activity relationships for cannabinoid receptor-binding and analgesic activity: Studies of bicyclic cannabinoid analogs. *Mol. Pharmacol.* **44**(5), 1008-1015 (1993).
2. Lindigkeit, R., Boehme, A., Eiserloh, I., *et al.* Spice: A never ending story? *Forensic Sci. Int.* **191**(1-3), 58-63 (2009).

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, 01/13/2026

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