

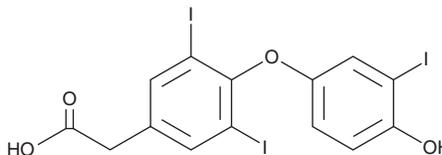
PRODUCT INFORMATION



Tiratricol

Item No. 28729

CAS Registry No.:	51-24-1
Formal Name:	4-(4-hydroxy-3-iodophenoxy)-3,5-diiodo-benzeneacetic acid
Synonym:	3,3',5-Triiodothyroacetic acid
MF:	C ₁₄ H ₉ I ₃ O ₄
FW:	621.9
Purity:	≥98%
UV/Vis.:	λ _{max} : 210, 227 nm
Supplied as:	A crystalline 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

Tiratricol is supplied as a crystalline solid. A stock solution may be made by dissolving the tiratricol in the solvent of choice, which should be purged with an inert gas. Tiratricol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of tiratricol in these solvents is approximately 30 mg/ml.

Description

Tiratricol is a metabolite of 3,3',5-triiodo-L-thyronine (Item No. 16028) and an agonist of the thyroid hormone receptors TR α and TR β (EC₅₀s = 1.81 and 4.13 nM, respectively, in reporter assays).¹ *In vivo*, tiratricol (50 or 400 ng/g) stimulates cerebellar Purkinje cell dendritogenesis and cortical myelination in *Mct8/Oatp1c1* double knockout mice.² It reduces goiter induced by methimazole (Item No. 23718) in rats when administered at a dose of 25 μ g per animal.³ Tiratricol also decreases toll-like receptor 2 (TLR2) levels and the phosphorylation of Akt and MAPK in HEK-Blue™ cells expressing human TLR2 and reduces liver tissue necrosis in a mouse model of hepatitis induced by concanavalin A (Item No. 14951).⁴

References

1. Oka, T., Mitsui-Watanabe, N., Tatarazako, N., *et al.* Establishment of transactivation assay systems using fish, amphibian, reptilian and human thyroid hormone receptors. *J. Appl. Toxicol.* **33(9)**, 991-1000 (2013).
2. Kersseboom, S., Horn, S., Visser, W.E., *et al.* In vitro and mouse studies supporting therapeutic utility of triiodothyroacetic acid in MCT8 deficiency. *Mol. Endocrinol.* **28(12)**, 1961-1970 (2014).
3. Alvarez, L., Burgueño, A., Zeni, S., *et al.* Comparison of the effects of 3,5,3'-triiodothyroacetic acid and triiodothyronine on goiter prevention and involution and on hepatic and skeletal parameters in rats. *Horm. Metab. Res.* **36(5)**, 291-297 (2004).
4. Ha, H.C., Jang, J.M., Zhou, D., *et al.* 3, 5, 3'-Triiodothyroacetic acid (TRIAC) is an anti-inflammatory drug that targets toll-like receptor 2. *Arch. Pharm. Res.* **41(10)**, 996-1008 (2018).

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, 12/19/2022

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