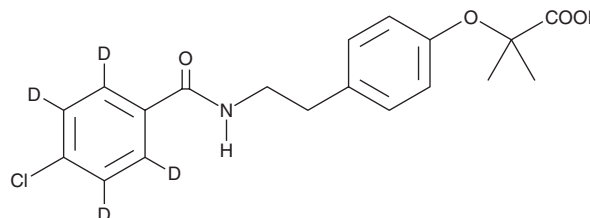


PRODUCT INFORMATION



Bezafibrate-d₄ Item No. 28526

CAS Registry No.: 1189452-53-6
Formal Name: 2-[4-[2-[(4-chlorobenzoyl-2,3,5,6-d₄)amino]ethyl]phenoxy]-2-methyl-propanoic acid
Synonym: Benzofibrate-d₄
MF: C₁₉H₁₆ClD₄NO₄
FW: 365.8
Chemical Purity: ≥98% (Bezafibrate)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
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

Bezafibrate-d₄ is intended for use as an internal standard for the quantification of bezafibrate (Item No. 10009145) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Bezafibrate-d₄ is supplied as a solid. A stock solution may be made by dissolving the bezafibrate-d₄ in the solvent of choice, which should be purged with an inert gas. Bezafibrate-d₄ is slightly soluble in DMSO.

Description

Bezafibrate is a non-selective agonist of peroxisome proliferator-activated receptors (PPARs; EC₅₀s = 50, 60, and 20 μM for human PPARα, PPARγ, and PPARδ, respectively).¹ It reduces triglyceride levels and the size of lipid droplets in an oleic acid-induced HepaRG hepatocyte model of steatosis when used at a concentration of 25 μM.² Bezafibrate (10 mg/kg per day) reduces plasma VLDL and LDL mass and triglyceride and free fatty acid levels in a high-fructose plus lard diet-induced rat model of insulin resistance.³

References

- Willson, T.M., Brown, P.J., Sternbach, D.D., *et al.* The PPARs: From orphan receptors to drug discovery. *J. Med. Chem.* **43**(4), 527-550 (2000).
- Rogue, A., Anthérieu, S., Vluggens, A., *et al.* PPAR agonists reduce steatosis in oleic acid-overloaded HepaRG cells. *Toxicol. Appl. Pharmacol.* **276**(1), 73-81 (2014).
- Matsui, H., Okumura, K., Kawakami, K., *et al.* Improved insulin sensitivity by bezafibrate in rats: Relationship to fatty acid composition of skeletal-muscle triglycerides. *Diabetes* **46**(3), 348-353 (1997).

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