

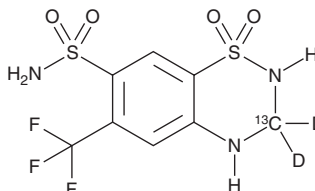
PRODUCT INFORMATION



Hydroflumethiazide-¹³C-d₂

Item No. 31608

CAS Registry No.: 1189877-11-9
Formal Name: 6-(trifluoromethyl)-3,4-dihydro-2H-benzo[e][1,2,4]thiadiazine-7-sulfonamide-3-¹³C-3,3-d₂ 1,1-dioxide
MF: C₇[¹³C]H₆D₂F₃N₃O₄S₂
FW: 334.3
Chemical Purity: ≥95% (Hydroflumethiazide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₂); ≤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

Hydroflumethiazide-¹³C-d₂ is intended for use as an internal standard for the quantification of hydroflumethiazide (Item No. 23612) 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).

Hydroflumethiazide-¹³C-d₂ is supplied as a solid. A stock solution may be made by dissolving the hydroflumethiazide-¹³C-d₂ in the solvent of choice, which should be purged with an inert gas. Hydroflumethiazide-¹³C-d₂ is soluble in methanol and DMSO.

Description

Hydroflumethiazide is a thiazide diuretic and an inhibitor of carbonic anhydrase (CA).¹ It is selective for CAII, CAVB, CAVII, CAIX, CAXII, and CAXIV (K_is = 0.24-0.31 μM) over CAI, CAIV, and CAVI (K_is = 2.8-8.3 μM), as well as CAIII, CAVA, and CAXIII (K_is = >10 μM). Formulations containing hydroflumethiazide have been used in the treatment of hypertension.

Reference

1. Temperini, C., Cecchi, A., Scozzafava, A., *et al.* Carbonic anhydrase inhibitors. Interaction of indapamide and related diuretics with 12 mammalian isozymes and X-ray crystallographic studies for the indapamide-isozyme II adduct. *Bioorg. Med. Chem. Lett.* **18**(8), 2567-2573 (2008).

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