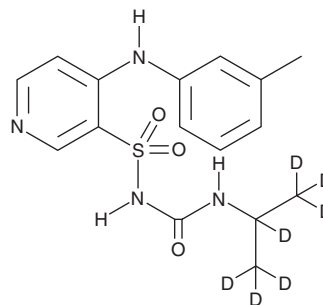


PRODUCT INFORMATION



Toraseamide-d₇ Item No. 26790

CAS Registry No.: 1189375-06-1
Formal Name: N-[[[1-methylethyl-d₇]amino]carbonyl]-4-[[3-methylphenyl]amino]-3-pyridinesulfonamide
Synonym: Torsemide-d₇
MF: C₁₆H₁₃D₇N₄O₃S
FW: 355.5
Chemical Purity: ≥98% (Toraseamide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤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

Toraseamide-d₇ is intended for use as an internal standard for the quantification of toraseamide 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).

Toraseamide-d₇ is supplied as a solid. A stock solution may be made by dissolving the toraseamide-d₇ in the solvent of choice, which should be purged with an inert gas. Toraseamide-d₇ is soluble in organic solvents such as methanol and DMSO.

Description

Toraseamide is a loop diuretic.¹ *In vivo*, toraseamide (6 mg/kg) increases fractional potassium, sodium, and chloride excretion and increases urine flow in rats. Formulations containing torsemide have been used to treat edema associated with congestive heart failure, kidney disease, and liver disease, as well as high blood pressure.

Reference

1. Beck, F.-X., Sone, M., Dorge, A., *et al.* Effect of loop diuretics on organic osmolytes and cell electrolytes in the renal outer medulla. *Kidney Int.* **42**(4), 843-850 (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, 12/13/2022

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