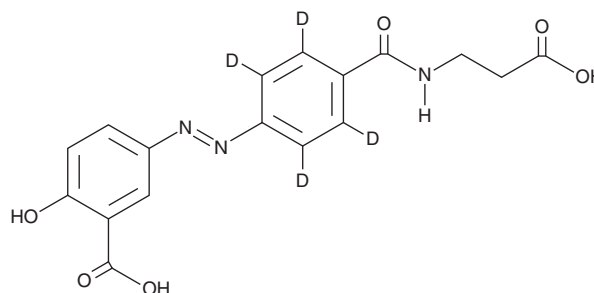


PRODUCT INFORMATION



Balsalazide-d₄ Item No. 28888

Formal Name: (E)-5-((4-((2-carboxyethyl)carbamoyl)phenyl-2,3,5,6-d₄)diazanyl)-2-hydroxybenzoic acid
MF: C₁₇H₁₁D₄N₃O₆
FW: 361.3
Chemical Purity: ≥98% (Balsalazide)
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

Balsalazide-d₄ is intended for use as an internal standard for the quantification of balsalazide (Item No. 18680) 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).

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

Description

Balsalazide is a prodrug form of 5-aminosalicylic acid (5-ASA; Item No. 70265).¹ It is cleaved by bacterial azoreductases in the intestinal lumen to release 5-ASA. Balsalazide (600 mg/kg per day) decreases IL-2 levels and increases IL-6 levels in the serum and colonic mucosa membrane, as well as decreases micro- and macroscopic colonic damage, in a rat model of colitis induced by 2,4-dinitrochlorobenzene (DNCB).² Formulations containing balsalazide have been used in the treatment of ulcerative colitis.

References

1. Tursi, A. Balsalazide in treating colonic diseases. *Expert Opin. Drug Metab. Toxicol.* **5**(12), 1555-1563 (2009).
2. Lu, Y.-X. and Chen, X.-M. Effects of balsalazide on experimental colitis induced by 2,4-dinitrochlorobenzene in rats. *Zhongguo Xinyao Yu Linchuang Zazhi* **26**(12), 901-904 (2007).

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, 12/06/2022

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