

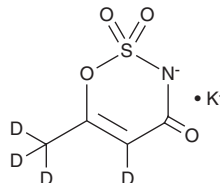
PRODUCT INFORMATION



Acesulfame-d₄ (potassium salt)

Item No. 27786

CAS Registry No.: 1623054-53-4
Formal Name: 6-(methyl-d₃)-1,2,3-oxathiazin-4(3H)-one-5-d,
2,2-dioxide, monopotassium salt
MF: C₄D₄NO₄S • K
FW: 205.3
Chemical Purity: ≥98% (Acesulfame)
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

Acesulfame-d₄ (potassium salt) is intended for use as an internal standard for the quantification of acesulfame 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).

Acesulfame-d₄ (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the acesulfame-d₄ (potassium salt) in the solvent of choice, which should be purged with an inert gas. Acesulfame-d₄ (potassium salt) is slightly soluble in methanol.

Description

Acesulfame is a non-caloric sweetener used as a sugar alternative that is not metabolized by humans.¹ Excreted acesulfame has been found in wastewater and is considered a pollutant.

Reference

1. Wu, M., Qian, Y., Boyd, J.M., *et al.* Direct large volume injection ultra-high performance liquid chromatography-tandem mass spectrometry determination of artificial sweeteners sucralose and acesulfame in well water. *J. Chromatogr. A.* **1359**, 156-161 (2014).

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, 11/02/2022

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