

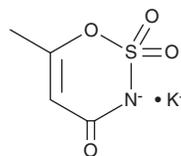
PRODUCT INFORMATION



Acesulfame (potassium salt)

Item No. 33356

CAS Registry No.: 55589-62-3
Formal Name: 6-methyl-1,2,3-oxathiazin-4(3H)-one
2,2-dioxide, monopotassium salt
Synonym: E950
MF: C₄H₄NO₄S • K
FW: 201.2
Purity: ≥98%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A crystalline 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 (potassium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the acesulfame (potassium salt) in the solvent of choice, which should be purged with an inert gas. Acesulfame (potassium salt) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of acesulfame (potassium salt) can be prepared by directly dissolving the acesulfame (potassium salt) in aqueous buffers. The solubility of acesulfame (potassium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

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, 11/01/2022

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