

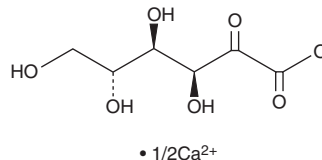
PRODUCT INFORMATION



2-keto D-Gluconic Acid (calcium salt)

Item No. 35622

CAS Registry No.: 3470-37-9
Formal Name: D-arabino-2-hexulosonic acid, hemicalcium salt
Synonyms: 2-keto D-Gluconate, 2-KGA
MF: $C_6H_9O_7 \cdot 1/2Ca$
FW: 213.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-keto D-Gluconic acid (calcium salt) is supplied as a solid. A stock solution may be made by dissolving the 2-keto D-gluconic acid (calcium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

2-keto D-Gluconic acid is a bacterial metabolite that has been found in *Pseudomonas* and a precursor in the synthesis of the antioxidant D-erythorbic acid.^{1,2}

References

1. Sun, L., Wang, D., Sun, W., *et al.* Characterization of a transcriptional regulator PtxS from *Pseudomonas plecoglossicida* for regulating 2-ketogluconic acid metabolism. *Int. J. Biol. Macromol.* **174**, 330-338 (2021).
2. Zeng, W., Cai, W., Liu, L., *et al.* Efficient biosynthesis of 2-keto-D-gluconic acid by fed-batch culture of metabolically engineered *Gluconobacter japonicus*. *Synth. Syst. Biotechnol.* **4**(3), 134-141 (2019).

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, 10/21/2022

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