

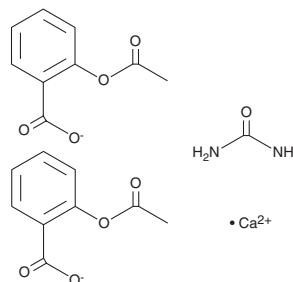
PRODUCT INFORMATION



Carbasalate Calcium

Item No. 29806

CAS Registry No.: 5749-67-7
Formal Name: 2-(acetyloxy)-benzoic acid, calcium salt, compd. with urea (2:1:1)
Synonym: Calcium Carbaspirin
MF: $C_{18}H_{14}O_8 \cdot Ca \cdot CH_4N_2O$
FW: 458.4
Purity: $\geq 95\%$
Supplied as: A crystalline 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

Carbasalate calcium is supplied as a crystalline solid. A stock solution may be made by dissolving the carbasalate calcium in the solvent of choice, which should be purged with an inert gas. Carbasalate calcium is slightly soluble ethanol, DMSO, and dimethyl formamide.

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 carbasalate calcium can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of carbasalate calcium in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Carbasalate calcium is a calcium-urea chelate of aspirin (Item No. 70260) that has analgesic, anti-inflammatory, and antipyretic effects.^{1,2} It reduces carrageenan-induced paw edema in rats.²

References

1. Murray, F.E., Hudson, N., Atherton, J.C., et al. Comparison of effects of calcium carbasalate and aspirin on gastroduodenal mucosal damage in human volunteers. *Gut* **38**(1), 11-14 (1996).
2. Liu, G., Han, Y., and Fan, X. Study on antipyretic, analgesic, anti-inflammatory and toxic effects of carbasalate calcium. *Zhongguo Yiyao Gongye Zazhi* **28**(10), 452-453 (1997).

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, 12/12/2024

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