

PRODUCT INFORMATION



Se-Aspirin

Item No. 19315

CAS Registry No.: 1850293-95-6
Formal Name: selenocyanic acid,
2-[[2-(acetyloxy)benzoyl]amino]
ethyl ester

Synonym: Selenium-Acetylsalicylic Acid

MF: C₁₂H₁₂N₂O₃Se

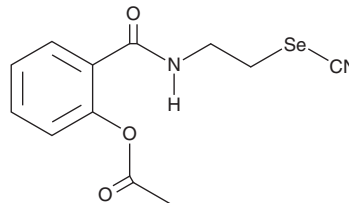
FW: 311.2

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

Se-Aspirin is supplied as a crystalline solid. A stock solution may be made by dissolving the Se-Aspirin in the solvent of choice. Se-Aspirin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of Se-Aspirin in these solvents is approximately 10, 30, and 50 mg/ml, respectively.

Se-Aspirin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Se-Aspirin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Se-Aspirin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Se-Aspirin is a hybrid of selenium and a nonsteroidal anti-inflammatory drug that has been shown to reduce the viability of different cancer cell lines, particularly colorectal cancer (CRC) cells (IC₅₀ = 3.4 μM).¹ It can inhibit the cell cycle in G₁ and G₂/M phases and induce apoptosis by activating caspase 3/7 and PARP cleavage.¹ Long-term exposure to Se-Aspirin is reported to cause an increase in intracellular reactive oxygen species levels in CRC cells.¹

Reference

1. Plano, D., Karelia, D.N., Pandey, M.K., *et al.* Design, synthesis, and biological evaluation of novel selenium (Se-NSAID) molecules as anticancer agents. *J. Med. Chem.* **59**(5), 1946-1959 (2016).

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, 06/06/2016

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