

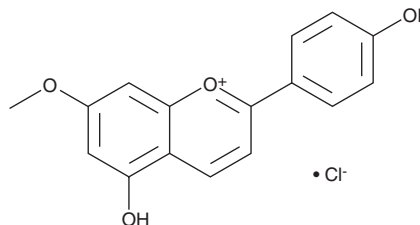
PRODUCT INFORMATION



7-methoxy Apigeninidin (chloride)

Item No. 19757

CAS Registry No.: 161773-51-9
Formal Name: 5-hydroxy-2-(4-hydroxyphenyl)-7-methoxy-1-benzopyrylium, monochloride
MF: C₁₆H₁₃O₄ • Cl
FW: 304.7
Purity: ≥98%
UV/Vis.: λ_{max}: 280, 481 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

7-methoxy Apigeninidin (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-methoxy apigeninidin (chloride) in the solvent of choice, which should be purged with an inert gas. 7-methoxy Apigeninidin (chloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 7-methoxy apigeninidin (chloride) in these solvents is approximately 1 mg/ml.

Description

7-methoxy Apigeninidin is a 3-deoxyanthocyanin fungicide originally isolated from *Sorghum* and a derivative of apigeninidin (Item No. 19756).^{1,2} It is active against the radial growth of *G. sorghi* and spore germination of *P. oryzae* (IC₅₀ = 50 ppm for both).²

References

1. Wu, X. and Prior, R.L. Identification and characterization of anthocyanins by high-performance liquid chromatography-electrospray ionization-tandem mass spectrometry in common foods in the United States: Vegetables, nuts, and grains. *J. Agric. Food Chem.* **53**(8), 3101-3113 (2005).
2. Aida, Y., Tamogami, S., Kodama, O., et al. Synthesis of 7-methoxyapigeninidin and its fungicidal activity against *Gloeocercospora sorghi*. *Biosci. Biotechnol. Biochem.* **60**(9), 1495-1496 (1996).

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, 12/19/2022

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