PRODUCT INFORMATION



Peonidin (chloride)

Item No. 19754

CAS Registry No.: Formal Name:	134-01-0 3,5,7-trihydroxy-2-(4-hydroxy-3- methoxyphenyl)-1-benzopyrylium, monochloride	ОН
MF:	C ₁₆ H ₁₃ O ₆ • Cl	HOOt
FW:	336.7	
UV/Vis.:	$\lambda_{\rm max}$: 276, 547 nm	OH •CI
Supplied as:	A crystalline solid	 ОН
Stability:	≥4 years	

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

Laboratory Procedures

Peonidin (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the peonidin (chloride) in the solvent of choice, which should be purged with an inert gas. Peonidin (chloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of peonidin (chloride) in ethanol and DMSO is approximately 16 mg/ml and approximately 25 mg/ml in DMF.

Peonidin (chloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, peonidin (chloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Peonidin (chloride) has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Peonidin is an O-methylated anthocyanidin that functions as a primary plant pigment, endowing purplish-red hues to flowers such as the peony, from which it takes its name, as well as berries and vegetables. It has been shown to exhibit chemopreventive, as well as anti-inflammatory activities on cancer cells in vitro, blocking COX-2 expression and transformation in JB6 P+ mouse epidermal cells.¹

Reference

1. Kwon, J. Y., Lee, K. W., Hur, H.J., et al. Peonidin inhibits phorbol-ester-induced COX-2 expression and transformation in JB6 P+ cells by blocking phosphorylation of ERK-1 and -2. Ann. N. Y. Acad. Sci. 1095, (2007).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 12/07/2022

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