

PRODUCT INFORMATION

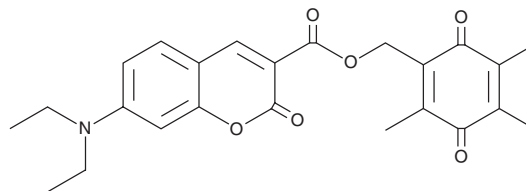


Coumarin-Quinone Conjugate

Item No. 29554

CAS Registry No.: 2741190-31-6
Formal Name: 7-(diethylamino)-2-oxo-2H-1-benzopyran-3-carboxylic acid, (2,4,5-trimethyl-3,6-dioxo-1,4-cyclohexadien-1-yl)methyl ester

MF: C₂₄H₂₅NO₆
FW: 423.5
Purity: ≥95%
UV/Vis.: λ_{max}: 212, 259, 421 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Coumarin-quinone conjugate is supplied as a solid. A stock solution may be made by dissolving the coumarin-quinone conjugate in the solvent of choice, which should be purged with an inert gas. Coumarin-quinone conjugate is soluble in DMSO and ethanol.

Description

Coumarin-quinone conjugate is a fluorescent substrate for NADH:ubiquinone oxidoreductases.¹ It is comprised of a coumarin fluorophore conjugated to a ubiquinone analog that can be reduced by NADH:ubiquinone oxidoreductases. It has been used to measure the kinetic parameters of apoptosis-inducing factor mitochondria-associated 2/ferroptosis suppressor protein 1 (AIFM2/FSP1).

Reference

1. Doll, S., Freitas, F.P., Shah, R., *et al.* FSP1 is a glutathione-independent ferroptosis suppressor. *Nature* **575(7784)**, 693-698 (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, 06/24/2025

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