

PRODUCT INFORMATION



DCI-Ac-Py

Item No. 39887

Formal Name: 2-bromo-4-((E)-2-(3-(dicyanomethylene)-5,5-dimethylcyclohex-1-en-1-yl)vinyl)-6-((E)-3-ethoxy-3-oxoprop-1-en-1-yl)phenyl isonicotinate

MF: $C_{30}H_{26}BrN_3O_4$

FW: 572.5

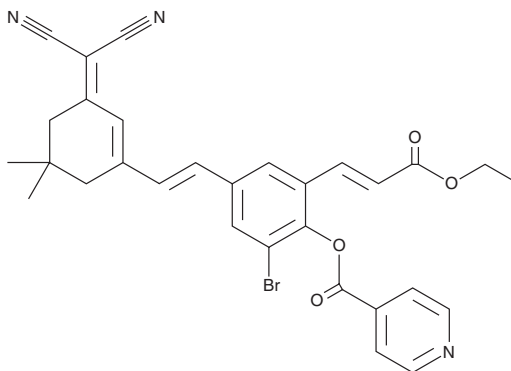
Purity: $\geq 98\%$

Abs./Em. MAX: 541/713 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

DCI-Ac-Py is supplied as a solid. A stock solution may be made by dissolving the DCI-Ac-Py in the solvent of choice, which should be purged with an inert gas. DCI-Ac-Py is soluble (≥ 10 mg/ml) in DMSO.

Description

DCI-Ac-Py is a turn-on fluorescent probe for the detection of biothiols.¹ Upon reaction with biothiols, it is converted into DCI-Br-Cou and displays a shift in absorbance from 396 to 541 nm and an emission maximum of 713 nm. DCI-Ac-Py has been used in the study of biothiol fluctuations in the brain in a mouse model of cerebral ischemia-reperfusion injury.

Reference

1. Yang, Y., Ma, M., Shen, L., *et al.* A fluorescent probe for investigating the role of biothiols in signaling pathways associated with cerebral ischemia-reperfusion injury. *Angew Chem. Int. Ed. Engl.* **62**, e2002310408 (2023).

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, 06/09/2025

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