

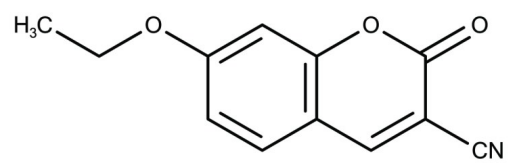
3-Cyano-7-ethoxycoumarin

Fluorescent substrate for Cytochrome P450

Fluorescent substrate for cytochrome P450. Excitation/emission = 408/455 nm. Can be used for CYP1A2 and CYP2C19 in intact cells. Cell permeable.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored unopened at room temperature
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Pale yellow solid.
CAS	117620-77-6
Couple Target	Cytochrome P450
Couple Type	Substrate
Emission Maximum	455 nm
Excitation Maximum	408 nm
Formula	$C_{12}H_9O_3$
Formulation	Lyophilized.
Identity	Identity determined by NMR.
MW	215.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (20 mg/ml, warm) or ethanol (3 mg/ml, warm).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com

Last modified: May 29, 2024