

PAP-AMC, Tyrosinase substrate

Fluorogenic substrate for tyrosinase. When assaying in phosphate buffer pH 7.1, the optimal excitation/emission wavelengths are 340/440 nm. 325/460 nm also work, but 380/460 nm will not detect this substrate.

Citations: 1

[View Online »](#)

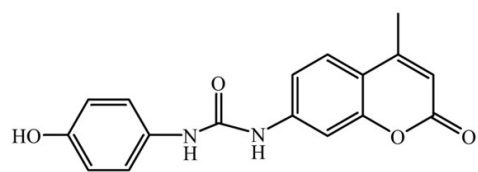
Ordering Information

[Order Online »](#)

BML-EI405-0100	0.1mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(4-Methyl-coumarin-7-yl)-3-(4-hydroxyphenyl)urea
Appearance	Off-white solid.
Couple Target	Tyrosinase
Couple Type	Substrate
Formula	$C_{17}H_{14}N_2O_4$
MW	310.3
MeltingPoint	238-242°C
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (18 mg/ml) or in methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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