

WP-AMC

DPP substrate

Fluorogenic substrate likely cleaved by dipeptidyl peptidase IV (DPIV, DPPIV, CD26), DPP8, and other prolyl peptidases. Can be used *in vitro*, for inhibitor screening, and in biological fluids. Ex = 380 nm, Em = 460 nm.

Ordering Information

[Order Online »](#)

BML-P450-0005	5mg
----------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Trp-Pro-AMC
Appearance	Off-white solid.
Formulation	Lyophilized.
Identity	Determined by MS.
MW	458.5
Purity	≥95% (HPLC)
Sequence	H-Trp-Pro-AMC
Solubility	Soluble in water or DMSO.

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)