

GP-AMC

DPP substrate

Fluorogenic substrate for dipeptidyl peptidase IV (DPIV, DPPIV, CD26) ($k_{cat}/K_m = 2.6 \times 10^5 \text{ s}^{-1} \text{ M}^{-1}$), DPPII/VII, DPP8 and DPP9, FAP (fibroblast activation protein), and likely POP/PREP, but not glutamate carboxypeptidase II. Can be used *in vitro*, for inhibitor screening, and in biological fluids. Ex = 380 nm, Em=460 nm.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P189-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Gly-Pro-AMC, Dipeptidyl peptidase substrate (fluorogenic)

CAS 115035-46-6

MW 410.3

Purity ≥97%

Sequence H-Gly-Pro-7-amino-4-methylcoumarin

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-
be@enzolifesciences.com](mailto:info-be@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)