

P32/98

DPPIV/DPP4 inhibitor

P32/98 (Ile-thiazolidide) is a specific, competitive transition-state substrate analog inhibitor of dipeptidyl peptidase IV (DPIV; DPPIV; CD26), with a K_i of 130 nM. Useful in diabetes research. Has been used *in vivo* and in tissue culture.

Citations: 5

[View Online »](#)

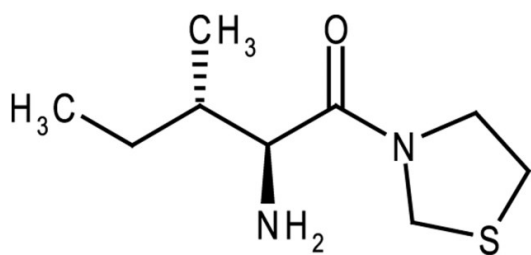
Ordering Information

[Order Online »](#)

BML-PI142-0010	10mg
BML-PI142-0050	50mg

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 3-N-(((2S,3S)-2-Amino-3-methylpentanoyl)-1,3-thiazolidine . hemi-fumarate

Appearance White solid.

Application Notes May be used to inhibit GLP-1 breakdown in blood, plasma and serum samples at a concentration of 1 mM.

CAS 136259-20-6

Couple Target DPP

Couple Type Inhibitor

Formula $C_9H_{18}N_2OS \cdot 0.5 C_4H_4O_4$

MW 260.4

Purity ≥98% (TLC, free base)

Solubility Soluble in water (20mg/ml) or DMSO (25mg/ml).

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)