

Diprotin A

DPP4 inhibitor

Inhibitor of dipeptidyl aminopeptidase IV.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-036-M025	25mg
------------------	------

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name H-Ile-Pro-Ile-OH

Appearance White to off-white solid.

Application Notes May be used to inhibit GLP-1 breakdown in blood, plasma and serum samples at a concentration of 100 µM.

CAS 90614-48-5

Formula $C_{17}H_{31}N_3O_4$

MW 341.5

Purity ≥98% (HPLC)

Solubility Soluble in water (1mg/ml).

Source Synthetic.

Last modified: May 29, 2024



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

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

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

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

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

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