

# 1G244

## DPP8/9 inhibitor

1G244 is a potent inhibitor of dipeptidyl peptidase 8 (DPP8) and dipeptidyl peptidase 9 (DPP9), with IC50 values of 14 nM and 53 nM, respectively.

Key features and applications include:

- **High Selectivity:** Specifically inhibits DPP8 and DPP9 without affecting DPPIV and DPPII.
- **Potency:** Demonstrates strong inhibitory effects with low IC50 values.
- **Induce Apoptosis:** Promotes apoptosis in multiple myeloma cells, exhibiting anti-myeloma effects.
- **Inhibit Adipogenesis:** Blocks adipogenesis in preadipocyte cell lines.
- **Suppress Tumor Growth:** Reduces subcutaneous tumor growth in murine xenograft models.

Relevant disease states include:

- **Multiple Myeloma:** Induces apoptosis in multiple myeloma cells, making it a valuable tool for studying this cancer.
- **Atherosclerosis:** Exhibits antiatherosclerotic properties by inhibiting DPP8/9.
- **Inflammatory Diseases:** Inhibition of DPP8/9 by 1G244 has been shown to cause pro-caspase-1-dependent pyroptosis in macrophages and monocytes.

## Ordering Information

[Order Online »](#)

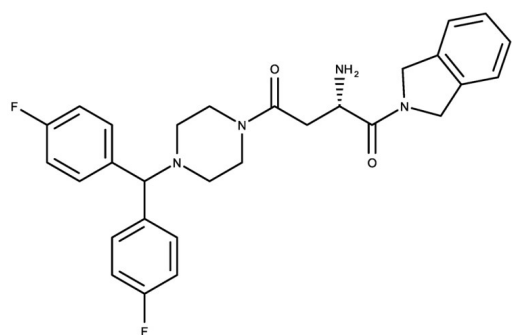
ENZ-CHM359-0025

25mg

## Manuals, SDS & CofA

[View Online »](#)

- GMP format available



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 2 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	(S)-2-Amino-4-{4-[bis-(4-fluorophenyl)-methyl]piperazin-1-yl}-1-(1,3-dihydro-isoindol-2-yl)-butane-1,4-dione, PTX-1210
Appearance	Off-white solid.
CAS	847928-32-9
Couple Target	DPP
Couple Type	Inhibitor
Formula	$C_{29}H_{30}F_2N_4O_2$
Identity	Determined by NMR.
MW	504.57
Purity	≥98% HPLC
Solubility	Soluble in DMSO (>20 mg/ml).



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)