

FABP4 inhibitor

A cell-permeable, potent, and selective inhibitor of FABP4.

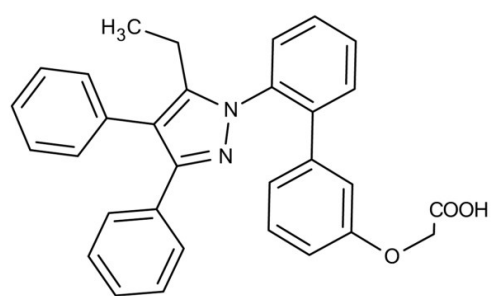
Ordering Information

[Order Online »](#)

ENZ-CHM118-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Adipocyte fatty-acid-binding protein inhibitor, ((2'-(5-Ethyl-3,4-diphenyl-1H-pyrazol-1-yl)(1,1'-biphenyl)-3-yl)oxy)-acetic acid, BMS 309403
Appearance	White powder solid.
CAS	300657-03-8
Couple Target	FABP
Couple Type	Inhibitor
Formula	$C_{31}H_{26}N_2O_3$
MW	474.6
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or ethanol.

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-fr@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