

# PF-622

## FAAH inhibitor

Novel, potent and selective inhibitor of fatty acid amide hydrolase (FAAH). Inhibits FAAH in a time-dependent manner by covalently modifying the enzyme's active site serine ( $IC_{50}=33nM$ ). It is completely selective for FAAH relative to other mammalian serine hydrolases.

Citations: 1

[View Online »](#)

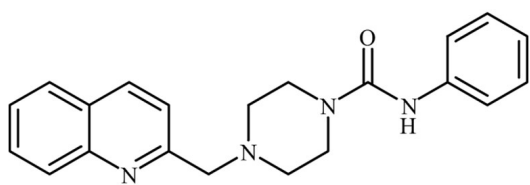
## Ordering Information

[Order Online »](#)

BML-EI402-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	N-phenyl-4-(quinolin-2-ylmethyl)piperazine-1-carboxamide
Appearance	Off-white solid.
CAS	898235-65-9
Couple Target	FAAH
Couple Type	Inhibitor
Formula	$C_{21}H_{22}N_4O$
MW	346.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (7mg/ml) or 100% ethanol (3mg/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-  
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)