

URB-602

Monoacylglycerol lipase inhibitor

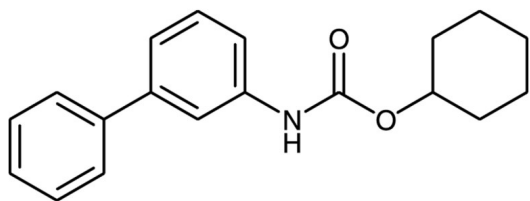
Selectively inhibits monoacylglycerol (MAG) lipase ($IC_{50}=28\mu M$) via a noncompetitive mechanism. Does not inhibit FAAH, DAG lipase or COX-2. URB-602 selectively raises levels of 2-arachidonoylglycerol with no change in anandamide concentrations in cells and tissues. In the same way anandamide levels were markedly elevated by the FAAH inhibitor URB597 at $1\mu M$.

Citations: 1

[View Online »](#)

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	Biphenyl-3-yl carbamic acid cyclohexyl ester
Appearance	Off-white solid.
CAS	565460-15-3
Couple Target	Monoacylglycerol lipase
Couple Type	Inhibitor
Formula	$C_{19}H_{21}NO_2$
MW	295.4
MeltingPoint	122-123°C
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (30 mg/ml) or ethanol (5 mg/ml warm).

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-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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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